

User Guide

# MILLIPLEX® Mouse Cytokine/Chemokine/ Growth Factor Expanded Magnetic Bead Panel 1

## 96-Well Plate Assay

**MCYT1-190K, MCYT1-190K-SPX, MCYT1-190K-SPXBK, MCYT1-190K-LPX, and  
MCYT1-190K-LPXBK**

|  |    |
|--|----|
| Introduction .....                                   | 3  |
| Principle .....                                      | 5  |
| Storage Conditions Upon Receipt .....                | 6  |
| Reagents Supplied.....                               | 7  |
| Mouse Cytokine/Chemokine/GF Expansion Panel 1        |    |
| Antibody-Immobilized Premixed Magnetic Beads .....   | 8  |
| Mouse Cytokine/Chemokine/GF Expansion Panel 1        |    |
| Antibody-Immobilized Individual Magnetic Beads ..... | 9  |
| Materials Required.....                              | 14 |
| Reagents .....                                       | 14 |
| Instrumentation/Materials.....                       | 14 |
| Safety Precautions .....                             | 15 |
| Symbol Definitions.....                              | 15 |
| Technical Guidelines .....                           | 18 |
| Sample Collection and Storage.....                   | 20 |
| Preparation of Serum Samples .....                   | 20 |
| Preparation of Plasma Samples.....                   | 20 |
| Preparation of Tissue Culture Supernatant.....       | 20 |

|   |    |
|---|----|
| Preparation of Reagents for Immunoassay .....                                 | 21 |
| Preparation of Antibody-Immobilized Beads.....                                | 21 |
| Preparation of Quality Controls .....   | 21 |
| Preparation of Wash Buffer .....  | 21 |
| Preparation of Serum Matrix .....   | 21 |
| Preparation of Mouse Cytokine/Chemokine/GF Expansion<br>Panel 1 Standard..... | 22 |
| Preparation of Standards .....  | 23 |
| Immunoassay Procedure .....   | 27 |
| Plate Washing.....  | 29 |
| Solid Plate.....  | 29 |
| Equipment Settings .....  | 30 |
| Quality Controls .....  | 33 |
| Assay Characteristics .....   | 33 |
| Cross-Reactivity .....  | 33 |
| Assay Sensitivities .....   | 33 |
| Precision .....   | 36 |
| Accuracy .....  | 39 |
| Troubleshooting .....   | 43 |
| Product Ordering.....   | 47 |
| Well Map.....   | 55 |
| Notice .....  | 56 |
| Technical Assistance .....  | 56 |
| Standard Warranty .....   | 56 |
| Contact Information.....  | 56 |
| Safety Data Sheets (SDS).....   | 56 |

---

## Introduction

Mouse models are fundamental in the field of biomedical research, serving as essential tools in preclinical and translational studies across various fields such as immunology, cancer, and infectious diseases. Cytokines, chemokines, and growth factors are key mediators of immune system functions capable of signaling through autocrine, paracrine, and endocrine mechanisms. Their pleiotropic immunomodulatory properties allow these biomolecules to react to diverse stimuli and regulate the immune response, either promoting or inhibiting inflammation.

The MILLIPLEX® portfolio offers the broadest selection of analytes across a wide range of disease states and species. Once the analytes of interest have been identified, you can rely on the quality that we build into each kit to produce results you can trust. In addition to the assay characteristics listed in the protocol, other performance criteria evaluated during the verification process include cross-reactivity, dilution linearity, kit stability, and sample behavior (for example detectability and stability).

Each MILLIPLEX® panel and kit includes:

- Quality controls (QCs) provided to qualify assay performance.
- Comparison of standard (calibrator) and QC lots to a reference lot to ensure lot-to-lot consistency.
- Optimized serum matrix to mimic native analyte environment.
- Detection antibody cocktails designed to yield consistent analyte profiles within panel.

In addition, each kit meets stringent manufacturing criteria to ensure batch-to-batch reproducibility. The MILLIPLEX® Mouse Cytokine/Chemokine/Growth Factor Expanded Panel 1 thus enables you to focus on the modulation of cytokine expression in mouse research models. Coupled with the Luminex® xMAP® platform in a magnetic bead format, you receive the advantage of ideal speed and sensitivity, allowing quantitative multiplex detection of dozens of analytes simultaneously, which can dramatically improve productivity.

The MILLIPLEX® Mouse Cytokine/Chemokine/Growth Factor Expanded Panel 1 is part of the most versatile system available for cytokine and chemokine research. From our single to multiplex biomarker solutions, we partner with you to design, develop, analytically verify and build the most comprehensive library available for protein detection and quantitation.

---

MILLIPLEX® kits offer you:

- The ability to select a 41-plex or 68-plex premixed option or
- The ability to choose any combination of analytes from our panel of 68 analytes to design a configurable kit that better meets your needs.
- A convenient “all-in-one” box format that gives you the assurance that you will have all the necessary reagents you need to run your assay.

The MILLIPLEX® Mouse Cytokine/Chemokine/Growth Factor Expanded Panel 1 is a 68-plex kit to be used for the simultaneous quantification of any or all of the following analytes in serum, plasma, or cell culture supernatant samples: BAFF, Betacellulin, CHI3L1, CXCL16, EGF, Eotaxin/CCL11, Erythropoietin, Exodus-2/CCL21/6Ckine, FGF-2, FLT3L, Fractalkine/CX3CL1, G-CSF, GDF-15/MIC-1, GM-CSF, Granzyme B, IFN $\alpha$ , IFN $\beta$ , IFN $\gamma$ , IL-10, IL-11, IL-12 (p40), IL-12 (p70), IL-13, IL-15, IL-16, IL-17A, IL-17F, IL-18, IL-1 $\alpha$ , IL-1 $\beta$ , IL-2, IL-20, IL-22, IL-28B/IFN $\lambda$ 3, IL-3, IL-31, IL-33/NF-HEV (mature), IL-4, IL-5, IL-6, IL-7, IL-9, IP-10/CXCL10, KC/GRO $\alpha$ /CXCL1, LIF, LIX/CXCL5, MCP-1/CCL2, MCP-2/CCL8, MCP-3/CCL7, MCP-5/CCL12, M-CSF, MDC/CCL22, MIG/CXCL9, MIP-1 $\alpha$ /CCL3, MIP-1 $\beta$ /CCL4, MIP-2/CXCL2, MIP-3 $\alpha$ /CCL20, MIP-3 $\beta$ /CCL19, RANTES/CCL5, sCD137/4-1BB/TNFRSF9, sFas/TNFRSF6, sFasL, sICAM-1, sRAGE, TARC/CCL17, TECK/CCL25, TNF $\alpha$ , and VEGF-A.

**For Research Use Only. Not for Use in Diagnostic Procedures.**

**Please read entire protocol before use.**

**It is important to use same assay incubation conditions throughout your study.**

## Principle

MILLIPLEX® kits are based on the Luminex® xMAP® technology — one of the fastest growing and most respected multiplex technologies offering applications throughout the life sciences and capable of performing a variety of bioassays including immunoassays on the surface of fluorescent-coded magnetic beads known as MagPlex® microspheres.

- Luminex® uses proprietary techniques to internally color-code microspheres with two fluorescent dyes. Through precise concentrations of these dyes, distinctly colored bead sets of 500 5.6 µm polystyrene microspheres or 80 6.45 µm magnetic microspheres can be created, each of which is coated with a specific capture antibody.
- After an analyte from a test sample is captured by the bead, a biotinylated detection antibody is introduced.
- The reaction mixture is then incubated with Streptavidin-PE conjugate, the reporter molecule, to complete the reaction on the surface of each microsphere.
- The following Luminex® instruments can be used to acquire and analyze data using two detection methods:
- The Luminex® analyzers Luminex® 200™, FLEXMAP 3D®, and xMAP® INTELLIFLEX are flow cytometry-based instruments that integrate key xMAP® detection components, such as lasers, optics, advanced fluidics and high-speed digital signal processors.
- The Luminex® analyzer, MAGPIX®, is a CCD-based instrument that integrates key xMAP® capture and detection components with the speed and efficiency of magnetic beads.
- Each individual microsphere is identified, and the result of its bioassay is quantified based on fluorescent reporter signals. We combine the streamlined data acquisition power of Luminex® xPONENT® acquisition software with sophisticated analysis capabilities of Belysa® Immunoassay Curve Fitting Software, integrating data acquisition and analysis seamlessly with all Luminex® instruments.
- xMAP® INTELLIFLEX runs on INTELLIFLEX software for instrument control, run setup and generating high quality data with flexible output options. Data can be exported in xPONENT® style CSV files for compatibility with many existing analytical applications, or in the new, customizable INTELLIFLEX file format. The INTELLIFLEX file format is intended for flexibility and simplicity, allowing the user to freely select which data points to include and to reduce the time to analysis.

The capability of adding multiple conjugated beads to each sample results in the ability to obtain multiple results from each sample. Open-architecture xMAP® technology enables multiplexing of many types of bioassays reducing time, labor and costs over traditional methods.

---

## Storage Conditions Upon Receipt

- Recommended storage for kit components is 2-8 °C.
- For long-term storage, freeze reconstituted standards and controls at  $\leq -20$  °C. Avoid multiple ( $> 2$ ) freeze/thaw cycles.
- DO NOT FREEZE Antibody-Immobilized Beads, Detection Antibody, and Streptavidin-Phycoerythrin.

## Reagents Supplied

| Reagents   | Volume               | Quantity    | Catalogue No.   |
|--|----------------------|-------------|---|
| Mouse Cytokine/Chemokine/GF Expansion Panel 1 Standard                 | 1 vial               | Lyophilized | MCYT1-8190-1 (for configurable 35-plex)<br>or<br>MCYT1-8190-2 (for 41-plex and 68-plex) |
| Mouse Cytokine/Chemokine/GF Expansion Panel 1 Quality Controls 1 and 2 | 1 vial each          | Lyophilized | MCYT1-6190-1 (for configurable 35-plex)<br>or<br>MCYT1-6190-2 (for 41-plex and 68-plex) |
| Serum Matrix   | 1 vial               | Lyophilized | MXMSM-MCYT1   |
| Set of one 96-Well Plate with 2 sealers                                | 1 plate<br>2 sealers | -           | -   |
| Assay Buffer   | 1 bottle             | 30 mL       | L-AB  |
| 10X Wash Buffer  | 1 bottle             | 60 mL       | L-WB  |
| Mouse Cytokine/Chemokine/GF Expansion Panel 1 Detection Antibodies     | 1 bottle             | 3.2 mL      | MCYT1-1190-1 (for configurable 35-plex)<br>or<br>MCYT1-1190-2 (for 41-plex and 68-plex) |
| Streptavidin-Phycoerythrin   | 1 bottle             | 3.2 mL      | L-SAPE20  |
| Mixing Bottle ( <i>not provided with premixed panel</i> )              | 1 bottle             | -           | -   |

For research use only. Not for use in diagnostic procedures.

## Mouse Cytokine/Chemokine/GF Expansion Panel 1 Antibody-Immobilized Premixed Magnetic Beads

| <b>Bead/Analyte Name</b> | <b>Volume</b> | <b>Quantity</b> | <b>Catalogue No.</b> |
|--------------------------|---------------|-----------------|----------------------|
| Premixed 41-plex Beads   | 3.5 mL        | 1 bottle        | MCYT1PMXS-MG         |
| Premixed 68-plex Beads   | 3.5 mL        | 1 bottle        | MCYT1PMXL-MG         |

Included Mouse Cytokine/Chemokine/GF Expansion Panel 1 Antibody-Immobilized Beads are dependent on customizable selection of analytes within the panel (see next page).



## Mouse Cytokine/Chemokine/GF Expansion Panel 1 Antibody-Immobilized Individual Magnetic Beads

**Configurable 68 Analytes  
(70X concentration, 70 µL)**

| <b>Bead/Analyte Name</b>        | <b>Luminex®<br/>Magnetic<br/>Bead Region</b> | <b>Instrument<br/>Compatibility<br/>Restrictions</b> | <b>41-Plex<br/>Magnetic<br/>Premixed<br/>Beads</b> | <b>68-Plex<br/>Magnetic<br/>Premixed<br/>Beads (not<br/>compatible<br/>with the<br/>MAGPIX®<br/>System)</b> | <b>Catalogue No.</b> |
|---------------------------------|--|--|--|---|----------------------|
| Anti-Mouse sFasL Bead           | 9  | Not compatible<br>with the<br>MAGPIX®<br>System      |  | ✓   | MSFASL-MG            |
| Anti-Mouse IL-22 Bead           | 12   | N/A  | ✓  | ✓   | MIL22-MG             |
| Anti-Mouse G-CSF Bead           | 13   | N/A  | ✓  | ✓   | MGCSF-MG             |
| Anti-Mouse<br>Eotaxin/CCL11Bead | 14   | N/A  | ✓  | ✓   | MCCL11-MG            |
| Anti-Mouse GM-CSF<br>Bead       | 15   | N/A  | ✓  | ✓   | MGMCSF-MG            |
| Anti-Mouse IL-33 Bead           | 18   | N/A  | ✓  | ✓   | MIL33-MG             |
| Anti-Mouse IFN $\gamma$ Bead    | 19   | N/A  | ✓  | ✓   | MIFNY-MG             |
| Anti-Mouse IL-1 $\alpha$ Bead   | 21   | N/A  | ✓  | ✓   | MIL1A-MG             |
| Anti-Mouse IL-17F Bead          | 22   | N/A  | ✓  | ✓   | MIL17F-MG            |
| Anti-Mouse IL-1 $\beta$ Bead    | 25   | N/A  | ✓  | ✓   | MIL1B-MG             |
| Anti- Mouse IL-2 Bead           | 26   | N/A  | ✓  | ✓   | MIL2-MG              |
| Anti-Mouse IFN $\beta$ Bead     | 27   | N/A  | ✓  | ✓   | MIFNB-MG             |
| Anti-Mouse IL-4 Bead            | 28   | N/A  | ✓  | ✓   | MIL4-MG              |
| Anti-Mouse IL-3 Bead            | 29   | N/A  | ✓  | ✓   | MIL3-MG              |
| Anti-Mouse IL-5 Bead            | 30   | N/A  | ✓  | ✓   | MIL5-MG              |
| Anti-Mouse IL-18 Bead           | 33   | N/A  | ✓  | ✓   | MIL18-MG             |
| Anti-Mouse IL-6 Bead            | 34   | N/A  | ✓  | ✓   | MIL6-MG              |

For research use only. Not for use in diagnostic procedures.

**Configurable 68 Analytes  
(70X concentration, 70 µL)**

| <b>Bead/Analyte Name</b>                  | <b>Luminex®<br/>Magnetic<br/>Bead Region</b> | <b>Instrument<br/>Compatibility<br/>Restrictions</b> | <b>41-Plex<br/>Magnetic<br/>Premixed<br/>Beads</b> | <b>68-Plex<br/>Magnetic<br/>Premixed<br/>Beads (not<br/>compatible<br/>with the<br/>MAGPIX®<br/>System)</b> | <b>Catalogue No.</b> |
|---|--|--|--|---|----------------------|
| Anti-Mouse FGF-2 Bead                     | 35   | N/A  | ✓  | ✓   | MFGF2-MG             |
| Anti-Mouse IL-7 Bead                      | 36   | N/A  | ✓  | ✓   | MIL7-MG              |
| Anti-Mouse EGF Bead                       | 37   | N/A  | ✓  | ✓   | MEGF-MG              |
| Anti-Mouse IL-9 Bead                      | 38   | N/A  | ✓  | ✓   | MIL9-MG              |
| Anti-Mouse IFN $\alpha$ Bead              | 39   | N/A  | ✓  | ✓   | MIFNA-MG             |
| Anti-Mouse<br>IL-28B/IFN $\lambda$ 3 Bead | 42   | N/A  | ✓  | ✓   | MIL28B-MG            |
| Anti-Mouse IL-10 Bead                     | 43   | N/A  | ✓  | ✓   | MIL10-MG             |
| Anti-Mouse MIP-3 $\alpha$<br>Bead         | 44   | N/A  |  | ✓   | MMIP3A-MG            |
| Anti-Mouse IL-12p40<br>Bead               | 45   | N/A  | ✓  | ✓   | MIL12P40-MG          |
| Anti-Mouse Granzyme<br>B Bead             | 46   | N/A  |  | ✓   | MGZMB-MG             |
| Anti-Mouse IL-12p70<br>Bead               | 47   | N/A  | ✓  | ✓   | MIL12P70-MG          |
| Anti-Mouse TARC Bead                      | 48   | N/A  |  | ✓   | MTARC-MG             |
| Anti-Mouse IL-16 Bead                     | 49   | Not compatible<br>with the<br>MAGPIX®<br>System      |  | ✓   | MIL16-MG             |
| Anti-Mouse LIF Bead                       | 51   | N/A  | ✓  | ✓   | MLIF-MG              |
| Anti-Mouse IL-13 Bead                     | 52   | N/A  | ✓  | ✓   | MIL13-MG             |
| Anti-Mouse LIX Bead                       | 53   | N/A  | ✓  | ✓   | MLIX-MG              |
| Anti-Mouse IL-15 Bead                     | 54   | N/A  | ✓  | ✓   | MIL15-MG             |

For research use only. Not for use in diagnostic procedures.

**Configurable 68 Analytes  
(70X concentration, 70 µL)**

| <b>Bead/Analyte Name</b>                  | <b>Luminex®<br/>Magnetic<br/>Bead Region</b> | <b>Instrument<br/>Compatibility<br/>Restrictions</b> | <b>41-Plex<br/>Magnetic<br/>Premixed<br/>Beads</b> | <b>68-Plex<br/>Magnetic<br/>Premixed<br/>Beads (not<br/>compatible<br/>with the<br/>MAGPIX®<br/>System)</b> | <b>Luminex®<br/>Magnetic<br/>Bead Region</b> |
|---|--|--|--|---|--|
| Anti-Mouse Exodus-2/<br>CCL21/6CKine Bead | 55   | N/A  |  | ✓   | MCCL21-MG                                    |
| Anti-Mouse IL-17A<br>Bead                 | 56   | N/A  | ✓  | ✓   | MIL17A-MG                                    |
| Anti-Mouse IP-10 Bead                     | 57   | N/A  | ✓  | ✓   | MIP10-MG                                     |
| Anti-Mouse Fas Bead                       | 58   | Not compatible<br>with the<br>MAGPIX®<br>System      |  | ✓   | MFAS-MG                                      |
| Anti-Mouse IL-20 Bead                     | 59   | Not compatible<br>with the<br>MAGPIX®<br>System      |  | ✓   | MIL20-MG                                     |
| Anti-Mouse KC Bead                        | 61   | N/A  | ✓  | ✓   | MKC-MG                                       |
| Anti-Mouse MCP-1 Bead                     | 62   | N/A  | ✓  | ✓   | MMCP1-MG                                     |
| Anti-Mouse MIP-3β<br>Bead                 | 63   | N/A  |  | ✓   | MMIP3B-MG                                    |
| Anti-Mouse MIP-1α<br>Bead                 | 64   | N/A  | ✓  | ✓   | MMIP1A-MG                                    |
| Anti-Mouse MDC Bead                       | 65   | N/A  |  | ✓   | MMDC-MG                                      |
| Anti-Mouse MIP-1β<br>Bead                 | 66   | N/A  | ✓  | ✓   | MMIP1B-MG                                    |
| Anti-Mouse M-CSF Bead                     | 67   | N/A  | ✓  | ✓   | MMCSF-MG                                     |
| Anti-Mouse sCD137/<br>4-1BB/TNFRSF9 Bead  | 68   | Not compatible<br>with the<br>MAGPIX®<br>System      |  | ✓   | MCD137-MG                                    |
| Anti-Mouse sICAM-1<br>Bead                | 69   | Not compatible<br>with the<br>MAGPIX®<br>System      |  | ✓   | MSICAM1-MG                                   |

For research use only. Not for use in diagnostic procedures.

**Configurable 68 Analytes  
(70X concentration, 70 µL)**

| <b>Bead/Analyte Name</b>   | <b>Luminex®<br/>Magnetic<br/>Bead Region</b> | <b>Instrument<br/>Compatibility<br/>Restrictions</b> | <b>41-Plex<br/>Magnetic<br/>Premixed<br/>Beads</b> | <b>68-Plex<br/>Magnetic<br/>Premixed<br/>Beads (not<br/>compatible<br/>with the<br/>MAGPIX®<br/>System)</b> | <b>Luminex®<br/>Magnetic<br/>Bead Region</b> |
|----------------------------|--|--|--|---|--|
| Anti-Mouse CXCL16 Bead     | 70   | Not compatible with the MAGPIX® System               |  | ✓   | MCXCL16-MG                                   |
| Anti-Mouse MCP-5 Bead      | 72   | N/A  |  | ✓   | MMCP5-MG                                     |
| Anti-Mouse MIP-2 Bead      | 73   | N/A  | ✓  | ✓   | MMIP2-MG                                     |
| Anti-Mouse MIG Bead        | 74   | N/A  | ✓  | ✓   | MMIG-MG                                      |
| Anti-Mouse RANTES Bead     | 75   | N/A  | ✓  | ✓   | MRANTES-MG                                   |
| Anti-Mouse VEGF-A Bead     | 76   | N/A  | ✓  | ✓   | MVEGFA-MG                                    |
| Anti-Mouse TNFα Bead       | 77   | N/A  | ✓  | ✓   | MTNFA-MG                                     |
| Anti-Mouse IL-11 Bead      | 78   | N/A  |  | ✓   | MIL11-MG                                     |
| Anti-Mouse sRAGE Bead      | 79   | Not compatible with the MAGPIX® System               |  | ✓   | MSRAGE-MG                                    |
| Anti-Mouse CCL25/TECK Bead | 80   | Not compatible with the MAGPIX® System               |  | ✓   | MCCL25-MG                                    |
| Anti-Mouse BAFF Bead       | 81   | Not compatible with the MAGPIX® System               |  | ✓   | MBAFF-MG                                     |
| Anti-Mouse GDF-15 Bead     | 82   | Not compatible with the MAGPIX® System               |  | ✓   | MGDF15-MG                                    |
| Anti-Mouse FLT3L Bead      | 83   | Not compatible with the MAGPIX® System               |  | ✓   | MFLT3L-MG                                    |

For research use only. Not for use in diagnostic procedures.

**Configurable 68 Analytes  
(70X concentration, 70 µL)**

| <b>Bead/Analyte Name</b>           | <b>Luminex®<br/>Magnetic<br/>Bead Region</b> | <b>Instrument<br/>Compatibility<br/>Restrictions</b> | <b>41-Plex<br/>Magnetic<br/>Premixed<br/>Beads</b> | <b>68-Plex<br/>Magnetic<br/>Premixed<br/>Beads (not<br/>compatible<br/>with the<br/>MAGPIX®<br/>System)</b> | <b>Luminex®<br/>Magnetic<br/>Bead Region</b> |
|------------------------------------|--|--|--|---|--|
| Anti-Mouse CHI3L1 Bead             | 84   | Not compatible with the MAGPIX® System               |  | ✓   | MCHI3L1-MG                                   |
| Anti-Mouse MCP-2 Bead              | 85   | Not compatible with the MAGPIX® System               |  | ✓   | MMCP2-MG                                     |
| Anti-Mouse MCP-3 Bead              | 86   | Not compatible with the MAGPIX® System               |  | ✓   | MMCP3-MG                                     |
| Anti-Mouse Erythropoietin Bead     | 87   | Not compatible with the MAGPIX® System               |  | ✓   | MEPO-MG                                      |
| Anti-Mouse Fractalkine/CX3CL1 Bead | 88   | Not compatible with the MAGPIX® System               |  | ✓   | MCX3CL1-MG                                   |
| Anti-Mouse IL-31 Bead              | 89   | Not compatible with the MAGPIX® System               |  | ✓   | MIL31-MG                                     |
| Anti-Mouse Betacellulin Bead       | 90   | Not compatible with the MAGPIX® System               |  | ✓   | MBTC-MG                                      |

For research use only. Not for use in diagnostic procedures.

---

## Materials Required (Not provided)

### Reagents

MAGPIX® Drive Fluid PLUS (Catalogue No. 40-50030), xMAP® Sheath Fluid PLUS (Catalogue No. 40-50021), or xMAP® Sheath Concentrate PLUS (Catalogue No. 40-50023).

### Instrumentation/Materials




- Adjustable Pipettes with Tips capable of delivering 25 µL to 1000 µL
- Multichannel Pipettes capable of delivering 5 µL to 50 µL, or 25 µL to 200 µL
- Reagent Reservoirs
- Polypropylene Microfuge Tubes
- Rubber Bands
- Aluminum Foil
- Absorbent Pads
- Laboratory Vortex Mixer
- Sonicator (Branson Ultrasonic Cleaner Model No. B200 or equivalent)
- Titer Plate Shaker (VWR® Microplate Shaker, Catalogue No. 12620-926 or equivalent)
- Luminex® 200™, FLEXMAP 3D®, MAGPIX® with xPONENT® software or xMAP® INTELLIFLEX with INTELLIFLEX software by Luminex® Corporation.
- Automatic Plate Washer for magnetic beads (BioTek® 405 LS and 405 TS, Catalogue. Nos. 40-094, 40-095, 40-096, 40-097 or equivalent) or Handheld Magnetic Separation Block (Catalogue No. 40-285 or equivalent).




**Note:** See Table on pages 8-11 for instrument requirements based on analyte selection.

## Safety Precautions







- All blood components and biological materials should be handled as potentially hazardous. Follow universal precautions as established by the Centers for Disease Control and Prevention and by the Occupational Safety and Health Administration when handling and disposing of infectious agents.
- Sodium azide or ProClin™ has been added to some reagents as a preservative. Although the concentrations are low, Sodium azide and ProClin™ may react with lead and copper plumbing to form highly explosive metal azides. Dispose of unused contents and waste in accordance with international, federal, state, and local regulations.

### Symbol Definitions

| Ingredient                                  | Catalogue No. | Label  |   |
|---|---------------|--|---|
| Mouse Cytokine Expansion Panel 1 Standard 1 | MCYT1-8190-1  |   |   |
| Mouse Cytokine Expansion Panel 1 Standard 2 | MCYT1-8190-2  | <br> | <p><b>Danger.</b> Harmful if swallowed, in contact with skin or if inhaled. Causes serious eye damage. May cause damage to organs Respiratory Tract through prolonged or repeated exposure. May cause damage to organs Brain through prolonged or repeated exposure if swallowed. Harmful to aquatic life with long lasting effects. Do not breathe dust. Wash skin thoroughly after handling. Do not eat, drink, or smoke when using this product. Use only outdoors or in a well-ventilated area. Avoid release to the environment. Wear protective gloves/ protective clothing/ eye protection/ face protection/ hearing protection. IF SWALLOWED: Call a POISON CENTER/ doctor if you feel unwell. Rinse mouth. IF ON SKIN: Wash with plenty of water. Call a POISON CENTER/ doctor if you feel unwell. IF INHALED: Remove person to fresh air and keep comfortable for breathing. Call a POISON CENTER/ doctor if you feel unwell. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. Get medical advice/ attention if you feel unwell. Take off contaminated clothing and wash it before reuse. Dispose of contents/ container to an approved waste disposal plant.</p> |

| Ingredient   | Catalogue No. | Label   |   |
|--|---------------|---|---|
| Mouse Cytokine Expansion Panel 1 QC 1 & 2 for Standard 1 | MCYT1-6190-1  |  |   |
| Mouse Cytokine Expansion Panel 1 QC 1 & 2 for Standard 2 | MCYT1-6190-2  |  |   |
|  |               |  | <p><b>Danger.</b> Harmful if swallowed, in contact with skin or if inhaled. Causes serious eye damage. May cause damage to organs Respiratory Tract through prolonged or repeated exposure. May cause damage to organs Brain through prolonged or repeated exposure if swallowed. Harmful to aquatic life with long lasting effects. Do not breathe dust. Wash skin thoroughly after handling. Do not eat, drink, or smoke when using this product. Use only outdoors or in a well-ventilated area. Avoid release to the environment. Wear protective gloves/ protective clothing/ eye protection/ face protection/ hearing protection. IF SWALLOWED: Call a POISON CENTER/ doctor if you feel unwell. Rinse mouth. IF ON SKIN: Wash with plenty of water. Call a POISON CENTER/ doctor if you feel unwell. IF INHALED: Remove person to fresh air and keep comfortable for breathing. Call a POISON CENTER/ doctor if you feel unwell. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. Get medical advice/ attention if you feel unwell. Take off contaminated clothing and wash it before reuse. Dispose of contents/ container to an approved waste disposal plant.</p> |
| Serum Matrix   | MXMSM-MCYT1   | No Label Required   | Harmful to aquatic life with long lasting effects. Avoid release to the environment. Dispose of contents/ container to an approved waste disposal plant.  |



| Ingredient  | Catalogue No. | Label  |  |
|---|---------------|--|--|
| Mouse Cytokine Expansion Panel 1 Detection Antibodies 1 | MCYT1-1190-1  |   | <b>Warning.</b> Causes serious eye irritation. May cause damage to organs Respiratory Tract through prolonged or repeated exposure. Do not breathe mist or vapours. Wash skin thoroughly after handling. Wear eye protection/ face protection. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. Get medical advice/ attention if you feel unwell. If eye irritation persists: Get medical advice/ attention. Dispose of contents/ container to an approved waste disposal plant. |
| Mouse Cytokine Expansion Panel 1 Detection Antibodies 2 | MCYT1-1190-2  |   | <b>Warning.</b> Causes serious eye irritation. May cause damage to organs Respiratory Tract through prolonged or repeated exposure. Do not breathe mist or vapours. Wash skin thoroughly after handling. Wear eye protection/ face protection. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. Get medical advice/ attention if you feel unwell. If eye irritation persists: Get medical advice/ attention. Dispose of contents/ container to an approved waste disposal plant. |
| Streptavidin-Phycoerythrin                              | L-SAPE20      | <br> | <b>Warning.</b> Causes serious eye irritation. May cause damage to organs Respiratory Tract through prolonged or repeated exposure. Do not breathe mist or vapours. Wash skin thoroughly after handling. Wear eye protection/ face protection. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing. Get medical advice/ attention if you feel unwell. If eye irritation persists: Get medical advice/ attention. Dispose of contents/ container to an approved waste disposal plant. |
| Assay Buffer  | L-AB          |   | <b>Warning.</b> Causes serious eye irritation. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.  |
| 10X Wash Buffer   | L-WB          |   | <b>Warning.</b> May cause an allergic skin reaction. Wear protective gloves. IF ON SKIN: Wash with plenty of soap and water.   |

## Technical Guidelines

To obtain reliable and reproducible results, the operator should carefully read this entire manual and fully understand all aspects of each assay step before running the assay. The following notes should be reviewed and understood before the assay is set up.

- FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.
- Do not use beyond the expiration date on the label.
- Do not mix or substitute reagents with those from other lots or sources.
- The Antibody-Immobilized Beads are light sensitive and must be protected from light at all times. Cover the assay plate containing beads with opaque plate lid or aluminum foil during all incubation steps.
- It is important to allow all reagents to warm to room temperature (20-25 °C) before use in the assay.
- Incomplete washing can adversely affect the assay outcome. All washing must be performed with the Wash Buffer provided.
- The standards prepared by serial dilution must be used within 1 hour of preparation. Discard any unused standards except the standard stock which may be stored at  $\leq -20$  °C for 1 month and at  $\leq -80$  °C for greater than one month.
- If samples fall outside the dynamic range of the assay, further dilute the samples with the appropriate diluent and repeat the assay.
- Any unused mixed Antibody-Immobilized Beads may be stored in the Mixing Bottle at 2-8 °C for up to one month.
- During the preparation of the standard curve, make certain to mix the higher concentration well before making the next dilution. Use a new tip with each dilution.
- The plate should be read immediately after the assay is finished. If, however, the plate cannot be read immediately, seal the plate, cover with aluminum foil or an opaque lid, and store the plate at 2-8 °C for up to 24 hours. Prior to reading, agitate the plate on the plate shaker at room temperature for 10 minutes. Delay in reading a plate may result in decreased sensitivity for some analytes.
- The titer plate shaker should be set at a speed to provide maximum orbital mixing without splashing of liquid outside the wells. For the recommended plate shaker, this would be a setting of 5-7 which is approximately 500-800 rpm.
- Ensure that the needle probe is clean. This may be achieved by sonication and/or alcohol flushes.
- When reading the assay on Luminex® 200™, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate or to the recommended filter plates using 3 alignment discs. When reading the assay on MAGPIX®, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate or to the recommended filter plates using 2 alignment discs. When reading the assay on FLEXMAP 3D®, adjust probe height according to the protocols recommended by Luminex® to the kit solid plate using 1 alignment disc.  
For FLEXMAP 3D® when using the solid plate in the kit, the final resuspension should be with 150  $\mu$ L Sheath Fluid PLUS in each well and 75  $\mu$ L should be aspirated.

For research use only. Not for use in diagnostic procedures.

---

For xMAP® INTELLIFLEX, adjust probe height based on the type of plate you are using, place an alignment disk or an alignment sphere in the well according to the protocol recommended by Luminex®.

- For cell culture supernatants or tissue extraction, use the culture or extraction medium as the matrix solution in background, standard curve and control wells. If samples are diluted in Assay Buffer, use the Assay Buffer as matrix.
- For serum/plasma samples that require further dilution beyond 1:2, use the serum matrix provided in the kit.
- For cell/tissue homogenate, the final cell or tissue homogenate should be prepared in a buffer that has a neutral pH, contains minimal detergents or strong denaturing detergents, and has an ionic strength close to physiological concentration. Avoid debris, lipids, and cell/tissue aggregates. Centrifuge samples before use.
- Vortex all reagents well before adding to plate.

# Sample Collection and Storage

## Preparation of Serum Samples

- Allow the blood to clot for at least 30 minutes before centrifugation for 10 minutes at 1000 x g. Remove serum and assay immediately or aliquot and store samples at  $\leq -20$  °C.
- Avoid multiple (> 2) freeze/thaw cycles.
- When using frozen samples, it is recommended to thaw the samples completely, mix well by vortexing and centrifuge prior to use in the assay to remove particulates.
- Serum samples should be diluted 1:2 in the Assay Buffer provided in the kit. For example, in a tube, 30  $\mu$ L of serum may be combined with 30  $\mu$ L of Assay Buffer. When further dilution beyond 1:2 is required, use Serum Matrix as the diluent.

## Preparation of Plasma Samples

- Plasma collection using EDTA as an anti-coagulant is recommended. Centrifuge for 10 minutes at 1000 x g within 30 minutes of blood collection. Remove plasma and assay immediately or aliquot and store samples at  $\leq -20$  °C.
- Avoid multiple (> 2) freeze/thaw cycles.
- When using frozen samples, it is recommended to thaw the samples completely, mix well by vortexing and centrifuge prior to use in the assay to remove particulates.
- Plasma samples should be diluted 1:2 in the Assay Buffer provided in the kit. For example, in a tube, 30  $\mu$ L of plasma may be combined with 30  $\mu$ L of Assay Buffer. When further dilution beyond 1:2 is required, use Serum Matrix as the diluent.

## Preparation of Tissue Culture Supernatant

Centrifuge the sample to remove debris and assay immediately or aliquot and store samples at  $\leq -20$  °C.

- Avoid multiple (> 2) freeze/thaw cycles.
- Tissue culture supernatant may require a dilution with an appropriate control medium prior to assay. Tissue/cell extracts should be done in neutral buffers containing reagents and conditions that do not interfere with assay performance. Excess concentrations of detergent, salt, denaturants, high or low pH, etc. will negatively affect the assay. Organic solvents should be avoided. The tissue/cell extract samples should be free of particles such as cells or tissue debris.

### **Note:**

- A maximum of 25  $\mu$ L per well of diluted serum or plasma can be used. Tissue culture or other media may also be used.
- All samples must be stored in polypropylene tubes. **DO NOT STORE SAMPLES IN GLASS.**
- Avoid debris, lipids and cells when using samples with gross hemolysis or lipemia.
- Care must be taken when using heparin as an anti-coagulant since an excess of heparin will provide falsely high values. Use no more than 10 IU heparin per mL of blood collected.

For research use only. Not for use in diagnostic procedures.

---

# Preparation of Reagents for Immunoassay

## Preparation of Antibody-Immobilized Beads

If **premixed beads** are used, sonicate the premixed bead bottle 30 seconds and then vortex for 1 minute before use.

For **individual vials of beads**, sonicate each antibody-bead vial for 30 seconds; vortex for 1 minute. Add 50  $\mu\text{L}$  from each antibody-bead vial to the Mixing Bottle and bring final volume to 3.5 mL with Assay Buffer. Vortex the mixed beads well. Unused portion may be stored at 2-8  $^{\circ}\text{C}$  for up to one month.

**Note:** Due to the composition of magnetic beads, you may notice a slight color in the bead solution. This does not affect the performance of the beads or the kit.

Example 1: When using 10 antibody-immobilized beads, add 50  $\mu\text{L}$  from each of the 10 bead vials to the Mixing Bottle. Then add 3.0 mL Assay Buffer.

Example 2: When using 35 antibody-immobilized beads, add 50  $\mu\text{L}$  from each of the 35 bead vials to the Mixing Bottle. Then add 1.75 mL Assay Buffer.

## Preparation of Quality Controls

Before use, reconstitute Quality Control 1 and Quality Control 2 with 250  $\mu\text{L}$  deionized water. Invert the vial several times to mix and vortex. Allow the vial to sit for 5-10 minutes. Transfer the reconstituted Quality Control 1 and Quality Control 2 into two polypropylene microfuge tubes. Unused portion may be stored at  $\leq -20^{\circ}\text{C}$  for up to one month.

## Preparation of Wash Buffer

Bring the 10X Wash Buffer to room temperature and mix to bring all salts into solution. Dilute 60 mL of 10X Wash Buffer with 540 mL deionized water. Store the unused portion at 2-8  $^{\circ}\text{C}$  for up to one month.

## Preparation of Serum Matrix

This step is required for serum or plasma samples only.

Add 2 mL Assay Buffer to the bottle containing lyophilized Serum Matrix. Mix well. Allow at least 10 minutes for complete reconstitution. Leftover reconstituted Serum Matrix should be stored at  $\leq -20^{\circ}\text{C}$  for up to one month.

## Preparation of Mouse Cytokine/Chemokine/GF Expansion Panel 1 Standard

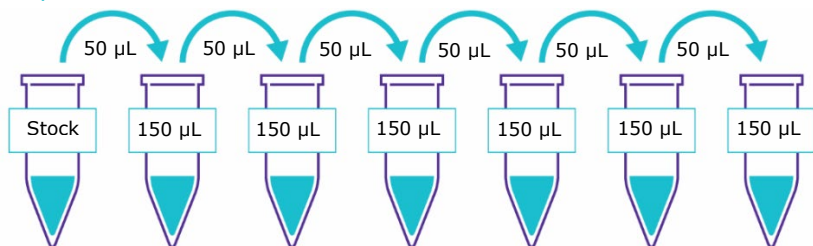
1. Prior to use, reconstitute the Mouse Cytokine/Chemokine/GF Expansion Panel 1 Standard with 250  $\mu\text{L}$  deionized water. Refer to table below for analyte concentrations. Invert the vial several times to mix. Vortex the vial for 10 seconds. Allow the vial to sit for 5-10 minutes. Transfer the reconstituted standard to a polypropylene microfuge tube. This will be used as Standard 7; the unused portion may be stored at  $\leq -20$   $^{\circ}\text{C}$  for up to one month.
2. Preparation of Working Standards

Label 6 polypropylene microfuge tubes Standard 1 through Standard 6. Add 150  $\mu\text{L}$  of Assay Buffer to each of the 6 tubes. Prepare serial dilutions by adding 50  $\mu\text{L}$  of the reconstituted standard to the Standard 6 tube, mix well and transfer 50  $\mu\text{L}$  of Standard 6 to the Standard 5 tube, mix well and transfer 50  $\mu\text{L}$  of Standard 5 to the Standard 4 tube, mix well and transfer 50  $\mu\text{L}$  of Standard 4 to the Standard 3 tube, mix well and transfer 50  $\mu\text{L}$  of Standard 3 to the Standard 2 tube, mix well and transfer 50  $\mu\text{L}$  of Standard 2 to the Standard 1 tube and mix well. The 0 Standard (Background) will be Assay Buffer.

| <b>Standard No.</b> | <b>Add Deionized Water (<math>\mu\text{L}</math>)</b> | <b>Add Standard (volume)</b> |
|---------------------|---|------------------------------|
| Standard 7          | 250 $\mu\text{L}$                                     | 0                            |

| <b>Standard No.</b> | <b>Add Assay Buffer (<math>\mu\text{L}</math>)</b> | <b>Add Standard (volume)</b>   |
|---------------------|--|--------------------------------|
| Standard 6          | 150 $\mu\text{L}$                                  | 50 $\mu\text{L}$ of Standard 7 |
| Standard 5          | 150 $\mu\text{L}$                                  | 50 $\mu\text{L}$ of Standard 6 |
| Standard 4          | 150 $\mu\text{L}$                                  | 50 $\mu\text{L}$ of Standard 5 |
| Standard 3          | 150 $\mu\text{L}$                                  | 50 $\mu\text{L}$ of Standard 4 |
| Standard 2          | 150 $\mu\text{L}$                                  | 50 $\mu\text{L}$ of Standard 3 |
| Standard 1          | 150 $\mu\text{L}$                                  | 50 $\mu\text{L}$ of Standard 2 |

## Preparation of Standards



Reconstituted Standard 7    Standard 6    Standard 5    Standard 4    Standard 3    Standard 2    Standard 1

| <b>Standard</b> | <b>IL-7 (pg/mL)</b> | <b>MIP-2, MCP-2 (pg/mL)</b> | <b>IL-3 (pg/mL)</b> |
|-----------------|---------------------|-----------------------------|---------------------|
| Standard 1      | 0.5                 | 0.6                         | 1.0                 |
| Standard 2      | 2.0                 | 2.4                         | 3.9                 |
| Standard 3      | 7.8                 | 10                          | 16                  |
| Standard 4      | 31                  | 39                          | 63                  |
| Standard 5      | 125                 | 156                         | 250                 |
| Standard 6      | 500                 | 625                         | 1,000               |
| Standard 7      | 2,000               | 2,500                       | 4,000               |

| <b>Standard</b> | <b>Fractalkine/CXC L1 (pg/mL)</b> | <b>GDF-15 (pg/mL)</b> | <b>MCP-3 (pg/mL)</b> |
|-----------------|-----------------------------------|-----------------------|----------------------|
| Standard 1      | 1.2                               | 1.5                   | 1.7                  |
| Standard 2      | 4.9                               | 5.9                   | 6.8                  |
| Standard 3      | 20                                | 23                    | 27                   |
| Standard 4      | 78                                | 94                    | 109                  |
| Standard 5      | 313                               | 375                   | 438                  |
| Standard 6      | 1,250                             | 1,500                 | 1,750                |
| Standard 7      | 5,000                             | 6,000                 | 7,000                |

For research use only. Not for use in diagnostic procedures.

| <b>Standard</b> | <b>IL-4, MDC<br/>(pg/mL)</b> | <b>GM-CSF (pg/mL)</b> | <b>IL-22, IL-17F,<br/>IL-2, IL-12p70,<br/>IL-17A,<br/>MIP-1<math>\alpha</math>,<br/>MCP-5,<br/>Betacellulin<br/>(pg/mL)</b> |
|-----------------|------------------------------|-----------------------|---|
|-----------------|------------------------------|-----------------------|---|

|            |       |       |        |
|------------|-------|-------|--------|
| Standard 1 | 1.8   | 2.2   | 2.4    |
| Standard 2 | 7.3   | 8.8   | 9.8    |
| Standard 3 | 29    | 35    | 39     |
| Standard 4 | 117   | 141   | 156    |
| Standard 5 | 469   | 563   | 625    |
| Standard 6 | 1,875 | 2,250 | 2,500  |
| Standard 7 | 7,500 | 9,000 | 10,000 |

| <b>Standard</b> | <b>Granzyme B<br/>(pg/mL)</b> | <b>IL-28B/IFN<math>\lambda</math>3,<br/>TARC (pg/mL)</b> | <b>Eotaxin/CCL11<br/>(pg/mL)</b> |
|-----------------|-------------------------------|--|----------------------------------|
|-----------------|-------------------------------|--|----------------------------------|

|            |        |        |        |
|------------|--------|--------|--------|
| Standard 1 | 2.9    | 3.1    | 3.2    |
| Standard 2 | 12     | 12     | 13     |
| Standard 3 | 47     | 49     | 51     |
| Standard 4 | 188    | 195    | 203    |
| Standard 5 | 750    | 781    | 813    |
| Standard 6 | 3,000  | 3,125  | 3,250  |
| Standard 7 | 12,000 | 12,500 | 13,000 |

| <b>Standard</b> | <b>IL-1<math>\alpha</math>, IL-5, IFN <math>\alpha</math>,<br/>MIP-3 <math>\alpha</math>, IL-13,<br/>VEGF-A, TNF <math>\alpha</math> (pg/mL)</b> | <b>IFN<math>\beta</math>, IL-9, IL-16,<br/>BAFF, FLT3L<br/>(pg/mL)</b> | <b>IFN<math>\gamma</math><br/>(pg/mL)</b> |
|-----------------|--|--|---|
|-----------------|--|--|---|

|            |        |        |        |
|------------|--------|--------|--------|
| Standard 1 | 3.7    | 4.9    | 6.1    |
| Standard 2 | 15     | 20     | 24     |
| Standard 3 | 59     | 78     | 98     |
| Standard 4 | 234    | 313    | 391    |
| Standard 5 | 938    | 1,250  | 1,563  |
| Standard 6 | 3,250  | 5,000  | 6,250  |
| Standard 7 | 15,000 | 20,000 | 25,000 |

For research use only. Not for use in diagnostic procedures.



| <b>Standard</b> | <b>IL-6, IL-12p40,<br/>Exodus-2/CCL21/<br/>6Ckine, RANTES<br/>(pg/mL)</b> | <b>LIX, IP-10<br/>(pg/mL)</b> | <b>LIF (pg/mL)</b> |
|-----------------|---|-------------------------------|--------------------|
| Standard 1      | 7.3   | 8.5                           | 9.8                |
| Standard 2      | 29  | 34                            | 39                 |
| Standard 3      | 117   | 137                           | 156                |
| Standard 4      | 469   | 547                           | 625                |
| Standard 5      | 1,875   | 2,188                         | 2,500              |
| Standard 6      | 7,500   | 8,750                         | 10,000             |
| Standard 7      | 30,000  | 35,000                        | 40,000             |

| <b>Standard</b> | <b>IL-1<math>\beta</math>, FGF-2,<br/>MCP-1, CXCL16,<br/>Erythropoietin<br/>(pg/mL)</b> | <b>IL-10, MIP-3<math>\beta</math>,<br/>sCD137/4-1BB/<br/>TNFRSF9, IL-11<br/>(pg/mL)</b> | <b>IL-18, Fas,<br/>IL-20, MIP-1<math>\beta</math>,<br/>M-CSF<br/>(pg/mL)</b> |
|-----------------|---|---|--|
| Standard 1      | 12  | 15  | 18   |
| Standard 2      | 49  | 59  | 73   |
| Standard 3      | 195   | 234   | 293  |
| Standard 4      | 781   | 938   | 1,172  |
| Standard 5      | 3,125   | 3,750   | 4,688  |
| Standard 6      | 12,500  | 15,000  | 18,750   |
| Standard 7      | 50,000  | 60,000  | 75,000   |

| <b>Standard</b> | <b>MIG (pg/mL)</b> | <b>KC (pg/mL)</b> | <b>IL-33, IL-15,<br/>CHI3L1 (pg/mL)</b> |
|-----------------|--------------------|-------------------|---|
| Standard 1      | 21                 | 22                | 25                                      |
| Standard 2      | 83                 | 88                | 98                                      |
| Standard 3      | 332                | 352               | 391                                     |
| Standard 4      | 1,328              | 1,406             | 1,563                                   |
| Standard 5      | 5,313              | 5,625             | 6,250                                   |
| Standard 6      | 21,250             | 22,500            | 25,000                                  |
| Standard 7      | 85,000             | 90,000            | 100,000                                 |

For research use only. Not for use in diagnostic procedures.

| <b>Standard</b> | <b>sFasL<br/>(pg/mL)</b> | <b>EGF, sRAGE,<br/>CCL25/TECK,<br/>IL-31 (pg/mL)</b> | <b>G-CSF<br/>(pg/mL)</b> | <b>sICAM-1<br/>(pg/mL)</b> |
|-----------------|--------------------------|--|--------------------------|----------------------------|
| Standard 1      | 37                       | 49   | 98                       | 122                        |
| Standard 2      | 147                      | 195  | 391                      | 488                        |
| Standard 3      | 586                      | 781  | 1,563                    | 1,953                      |
| Standard 4      | 2,344                    | 3,125  | 6,250                    | 7,813                      |
| Standard 5      | 9,375                    | 12,500   | 25,000                   | 31,250                     |
| Standard 6      | 37,500                   | 50,000   | 100,000                  | 125,000                    |
| Standard 7      | 150,000                  | 200,000  | 400,000                  | 500,000                    |

For research use only. Not for use in diagnostic procedures.

# Immunoassay Procedure

- Prior to beginning this assay, it is imperative to read this protocol completely and to thoroughly understand the Technical Guidelines.
- Allow all reagents to warm to room temperature (20-25 °C) before use in the assay.
- Diagram the placement of Standards, 0 (Background), Standard 1 through 7, Controls 1 and 2, and Samples on Well Map Worksheet in a vertical configuration. It is recommended to run the assay in duplicate.

**Note:** Most instruments will only read the 96-well plate vertically by default.

1. Add 200  $\mu\text{L}$  of Wash Buffer into each well of the plate. Seal and mix on a plate shaker for 10 minutes at room temperature (20-25 °C).
2. Decant Wash Buffer and remove the residual amount from all wells by inverting the plate and tapping it smartly onto absorbent towels several times.
3. Add 25  $\mu\text{L}$  of each Standard or Control into the appropriate wells. Assay Buffer should be used for 0 standard (Background).
4. Add 25  $\mu\text{L}$  of Assay Buffer to the sample wells.
5. Add 25  $\mu\text{L}$  of appropriate matrix solution to the background, standards, and control wells. When assaying serum or plasma, use the Serum Matrix. When assaying tissue culture or other supernatant, use proper control culture medium as the matrix solution.
6. Add 25  $\mu\text{L}$  of Sample (1:2 diluted or neat) into the appropriate wells.
7. Vortex Mixing Bottle and add 25  $\mu\text{L}$  of the Mixed or Premixed Beads to each well.

**Note:** During addition of Beads, shake bead bottle intermittently to avoid settling.

8. Seal the plate with a plate sealer. Wrap the plate with foil and incubate with agitation on a plate shaker overnight (16-18 hours) at 2-8 °C. Alternatively, incubate for 2 hours at room temperature (20-25 °C).

Add 200  $\mu\text{L}$  Wash Buffer per well



Shake 10 min, RT  
Decant

- Add 25  $\mu\text{L}$  Standard or Control to appropriate wells
- Add 25  $\mu\text{L}$  Assay Buffer to background and sample wells
- Add 25  $\mu\text{L}$  appropriate matrix solution to background, standards, and control wells
- Add 25  $\mu\text{L}$  1:2 diluted Samples to sample wells
- Add 25  $\mu\text{L}$  Beads to each well



Incubate overnight (16-18 hours) at 2-8 °C or 2 hours at 20-25 °C.

9. Gently remove well contents and wash plate 3 times following instructions listed in the PLATE WASHING section.
10. Add 25  $\mu\text{L}$  of Detection Antibodies into each well.  
**Note:** Allow the Detection Antibodies to warm to room temperature prior to addition.
11. Seal, cover with foil and incubate with agitation on a plate shaker for 1 hour at room temperature (20-25  $^{\circ}\text{C}$ ).
12. Gently remove well contents and wash plate 3 times following instructions listed in the PLATE WASHING section.
13. Add 25  $\mu\text{L}$  Streptavidin-Phycoerythrin to each well.
14. Seal, cover with foil and incubate with agitation on a plate shaker for 30 minutes at room temperature (20-25  $^{\circ}\text{C}$ ).
15. Gently remove well contents and wash plate 3 times following instructions listed in the PLATE WASHING section.
16. Add 150  $\mu\text{L}$  of Sheath Fluid PLUS (or Drive Fluid PLUS if using MAGPIX<sup>®</sup>) to all wells. Resuspend the beads on a plate shaker for 5 minutes.
17. Run plate on Luminex<sup>®</sup> 200<sup>™</sup>, HTS, FLEXMAP 3D<sup>®</sup>, MAGPIX<sup>®</sup> with xPONENT<sup>®</sup> software or xMAP<sup>®</sup> INTELLIFLEX with INTELLIFLEX Software.
18. Save and analyze the Median Fluorescent Intensity (MFI) data using a 5-parameter logistic or spline curve-fitting method for calculating analyte concentrations in samples.

**Note:** For diluted samples, final sample concentrations should be multiplied by the dilution factor. For samples diluted as per protocol instructions, multiply by 2. If using another dilution factor, multiply by the appropriate dilution factor.



Remove well contents and wash 3X with 200  $\mu\text{L}$  Wash Buffer

Add 25  $\mu\text{L}$  Detection Antibodies per well



Incubate 1 hour at RT  
Remove well contents and wash 3X with 200  $\mu\text{L}$  Wash Buffer

Add 25  $\mu\text{L}$  Streptavidin-Phycoerythrin per well



Incubate 30 minutes at RT

Remove well contents and wash 3X with 200  $\mu\text{L}$  Wash Buffer

Add 150  $\mu\text{L}$  Sheath Fluid PLUS or Drive Fluid PLUS per well

Read on Luminex<sup>®</sup> instrument (100  $\mu\text{L}$ , 50 beads per bead set)

---

## Plate Washing

If using a solid plate, use either a handheld magnet or magnetic plate washer.

### Solid Plate

- Handheld magnet (Catalogue No. 40-285)

Rest plate on magnet for 60 seconds to allow complete settling of magnetic beads. Remove well contents by gently decanting the plate in an appropriate waste receptacle and gently tapping on absorbent pads to remove residual liquid. Wash plate with 200  $\mu\text{L}$  of Wash Buffer by removing plate from magnet, adding Wash Buffer, shaking for 30 seconds, reattaching to magnet, letting beads settle for 60 seconds and removing well contents as previously described after each wash. Repeat wash steps as recommended in Assay Procedure.

- Magnetic plate washer (Catalogue Nos. 40-094, 40-095, 40-096 and 40-097)

Please refer to specific automatic plate washer manual for appropriate equipment settings. Please note that after the final aspiration, there will be approximately 25  $\mu\text{L}$  of residual wash buffer in each well. This is expected when using the BioTek® plate washer and this volume does not need to be aspirated from the plate.

If using an automatic plate washer other than BioTek® 405 LS or 405 TS, please refer to the manufacturer's recommendations for programming instructions.

## Equipment Settings

Luminex® 200™, FLEXMAP 3D®, MAGPIX® with xPONENT® software and xMAP® INTELLIFLEX with INTELLIFLEX software:

These specifications are for the above listed instruments and software. Luminex® instruments with other software (for example MasterPlex®, StarStation, LiquiChip, Bio-Plex® Manager™, LABScan™100) would need to follow instrument instructions for gate settings and additional specifications from the vendors for reading Luminex® magnetic beads.

For magnetic bead assays, each instrument must be calibrated, and performance verified with the indicated calibration and verification kits.

| Instrument        | Calibration Kit  | Verification Kit  |
|-------------------|--|---|
| Luminex® 200™     | xPONENT® 3.1 compatible Calibration Kit (Catalogue No. LX2R-CAL-K25) | Performance Verification Kit (Catalogue No. LX2R-PVER-K25)                  |
| FLEXMAP 3D®       | FLEXMAP 3D® Calibrator Kit (Catalogue No. F3D-CAL-K25)               | FLEXMAP 3D® Performance Verification Kit (Catalogue No. F3D-PVER-K25)       |
| xMAP® INTELLIFLEX | xMAP® INTELLIFLEX Calibration Kit (Catalogue No. IFX-CAL-K20)        | xMAP® INTELLIFLEX Performance Verification Kit (Catalogue No. IFX-PVER-K20) |
| MAGPIX®           | MAGPIX® Calibration Kit (Catalogue No. MPX-CAL-K25)                  | MAGPIX® Performance Verification Kit (Catalogue No. MPX-PVER-K25)           |

**Note:** When setting up a Protocol using the xPONENT® software, you must select MagPlex® as the Bead Type in the Acquisition settings.

**Note:** These assays cannot be run on any instruments using Luminex® IS 2.3 or Luminex® 1.7 software.

The Luminex® probe height must be adjusted to the plate provided in the kit. Please use Catalogue No. MAG-PLATE, if additional plates are required for this purpose.

|                        |                            |
|------------------------|----------------------------|
| Events                 | 50, per bead               |
| Sample Size            | 100 $\mu$ L                |
| Gate Settings          | 8,000 to 15,000            |
| Reporter Gain          | Default (low PMT)          |
| Time Out               | 60 seconds                 |
| Bead Set               | Customizable 68-plex Beads |
|                        | <hr/>                      |
| sFasL                  | 9                          |
| IL-22                  | 12                         |
| G-CSF                  | 13                         |
| Eotaxin/CCL11          | 14                         |
| GM-CSF                 | 15                         |
| IL-33                  | 18                         |
| IFN $\gamma$           | 19                         |
| IL-1 $\alpha$          | 21                         |
| IL-17F                 | 22                         |
| IL-1 $\beta$           | 25                         |
| IL-2                   | 26                         |
| IFN $\beta$            | 27                         |
| IL-4                   | 28                         |
| IL-3                   | 29                         |
| IL-5                   | 30                         |
| IL-18                  | 33                         |
| IL-6                   | 34                         |
| FGF-2                  | 35                         |
| IL-7                   | 36                         |
| EGF                    | 37                         |
| IL-9                   | 38                         |
| IFN $\alpha$           | 39                         |
| IL-28B/IFN $\lambda$ 3 | 42                         |

For research use only. Not for use in diagnostic procedures.

---

|                           |    |
|---------------------------|----|
| IL-10                     | 43 |
| MIP-3 $\alpha$            | 44 |
| IL-12p40                  | 45 |
| Granzyme B                | 46 |
| IL-12p70                  | 47 |
| TARC                      | 48 |
| IL-16                     | 49 |
| LIF                       | 51 |
| IL-13                     | 52 |
| LIX                       | 53 |
| IL-15                     | 54 |
| Exodus-2/<br>CCL21/6Ckine | 55 |
| IL-17A                    | 56 |
| IP-10                     | 57 |
| Fas                       | 58 |
| IL-20                     | 59 |
| KC                        | 61 |
| MCP-1                     | 62 |
| MIP-3 $\beta$             | 63 |
| MIP-1 $\alpha$            | 64 |
| MDC                       | 65 |
| MIP-1 $\beta$             | 66 |
| M-CSF                     | 67 |
| sCD137/4-1BB/<br>TNFRSF9  | 68 |
| sICAM-1                   | 69 |
| CXCL16                    | 70 |
| MCP-5                     | 72 |
| MIP-2                     | 73 |
| MIG                       | 74 |
| RANTES                    | 75 |



|                    |    |
|--------------------|----|
| VEGF-A             | 76 |
| TNF $\alpha$       | 77 |
| IL-11              | 78 |
| sRAGE              | 79 |
| CCL25/TECK         | 80 |
| BAFF               | 81 |
| GDF-15             | 82 |
| FLT3L              | 83 |
| CHI3L1             | 84 |
| MCP-2              | 85 |
| MCP-3              | 86 |
| Erythropoietin     | 87 |
| Fractalkine/CX3CL1 | 88 |
| IL-31              | 89 |
| Betacellulin       | 90 |

## Quality Controls

The ranges for each analyte in Quality Control 1 and 2 are provided on the card insert or can be located at our website [SigmaAldrich.com](https://www.sigmaaldrich.com) using the catalogue number as the keyword.

## Assay Characteristics

### Cross-Reactivity

There was no or negligible cross-reactivity between the antibodies for an analyte and any of the other analytes in this panel.

### Assay Sensitivities (minimum detectable concentrations, pg/mL)

Minimum Detectable Concentration (MinDC) is calculated using MILLIPLEX® Analyst 5.1. It measures the true limits of detection for an assay by mathematically determining what the empirical MinDC would be if an infinite number of standard concentrations were run for the assay under the same conditions.

| Analyte                | Overnight Protocol<br>(n = 13 Assays) |                      | 2-Hour Protocol<br>(n = 3 Assays) |                      |
|------------------------|---------------------------------------|----------------------|-----------------------------------|----------------------|
|                        | MinDC<br>(pg/mL)                      | MinDC+2SD<br>(pg/mL) | MinDC<br>(pg/mL)                  | MinDC+2SD<br>(pg/mL) |
| sFasL                  | 20.08                                 | 29.18                | 16.87                             | 32.02                |
| IL-22                  | 1.98                                  | 2.63                 | 1.36                              | 2.07                 |
| G-CSF                  | 43.62                                 | 75.60                | 66.67                             | 174.97               |
| Eotaxin/CCL11          | 2.75                                  | 3.61                 | 2.95                              | 3.38                 |
| GM-CSF                 | 0.58                                  | 0.84                 | 0.71                              | 1.18                 |
| IL-33                  | 11.09                                 | 19.25                | 18.07                             | 43.68                |
| IFN $\gamma$           | 1.46                                  | 2.08                 | 1.91                              | 3.08                 |
| IL-1 $\alpha$          | 1.37                                  | 2.31                 | 2.46                              | 3.23                 |
| IL-17F                 | 0.83                                  | 1.49                 | 1.35                              | 2.83                 |
| IL-1 $\beta$           | 6.61                                  | 10.66                | 4.49                              | 7.74                 |
| IL-2                   | 1.19                                  | 2.12                 | 0.98                              | 1.72                 |
| IFN $\beta$            | 2.52                                  | 4.03                 | 3.39                              | 5.36                 |
| IL-4                   | 0.52                                  | 0.74                 | 0.42                              | 0.54                 |
| IL-3                   | 0.28                                  | 0.51                 | 0.47                              | 0.84                 |
| IL-5                   | 2.07                                  | 4.04                 | 2.15                              | 4.82                 |
| IL-18                  | 16.15                                 | 23.60                | 9.10                              | 18.10                |
| IL-6                   | 2.49                                  | 4.17                 | 2.85                              | 6.68                 |
| FGF-2                  | 18.76                                 | 30.75                | 12.75                             | 37.73                |
| IL-7                   | 0.13                                  | 0.19                 | 0.21                              | 0.26                 |
| EGF                    | 28.51                                 | 43.53                | 27.16                             | 61.78                |
| IL-9                   | 3.16                                  | 6.10                 | 3.63                              | 6.81                 |
| IFN $\alpha$           | 2.09                                  | 3.52                 | 1.33                              | 2.56                 |
| IL-28B/IFN $\lambda$ 3 | 1.61                                  | 3.71                 | 2.56                              | 7.15                 |
| IL-10                  | 4.88                                  | 8.62                 | 6.87                              | 17.67                |
| MIP-3 $\alpha$         | 3.04                                  | 4.97                 | 3.15                              | 5.46                 |
| IL-12p40               | 2.57                                  | 4.91                 | 2.58                              | 3.80                 |
| Granzyme B             | 1.08                                  | 1.91                 | 1.41                              | 3.91                 |
| IL-12p70               | 0.86                                  | 1.44                 | 2.31                              | 5.58                 |

For research use only. Not for use in diagnostic procedures.

| <b>Analyte</b>            | <b>Overnight Protocol<br/>(n = 13 Assays)</b> |                              | <b>2-Hour Protocol<br/>(n = 3 Assays)</b> |                              |
|---------------------------|---|------------------------------|---|------------------------------|
|                           | <b>MinDC<br/>(pg/mL)</b>                      | <b>MinDC+2SD<br/>(pg/mL)</b> | <b>MinDC<br/>(pg/mL)</b>                  | <b>MinDC+2SD<br/>(pg/mL)</b> |
| TARC                      | 1.59  | 2.33                         | 2.07                                      | 4.04                         |
| IL-16                     | 4.02  | 5.75                         | 3.59                                      | 4.11                         |
| LIF                       | 6.29  | 9.57                         | 6.05                                      | 13.37                        |
| IL-13                     | 1.52  | 2.75                         | 2.47                                      | 3.94                         |
| LIX                       | 6.63  | 9.81                         | 6.19                                      | 6.66                         |
| IL-15                     | 12.26   | 22.18                        | 25.68                                     | 43.72                        |
| Exodus-2/<br>CCL21/6Ckine | 4.13  | 6.81                         | 5.70                                      | 6.67                         |
| IL-17A                    | 0.83  | 1.40                         | 1.03                                      | 2.46                         |
| IP-10                     | 6.60  | 8.26                         | 6.19                                      | 9.56                         |
| Fas                       | 7.44  | 12.83                        | 8.71                                      | 22.40                        |
| IL-20                     | 16.75   | 23.30                        | 10.61                                     | 29.30                        |
| KC                        | 12.18   | 26.87                        | 14.49                                     | 29.97                        |
| MCP-1                     | 4.98  | 9.83                         | 5.88                                      | 15.25                        |
| MIP-3 $\beta$             | 14.86   | 16.16                        | 13.96                                     | 16.30                        |
| MIP-1 $\alpha$            | 1.62  | 2.90                         | 1.44                                      | 3.06                         |
| MDC                       | 0.97  | 2.14                         | 0.87                                      | 1.70                         |
| MIP-1 $\beta$             | 17.28   | 24.90                        | 15.58                                     | 22.54                        |
| M-CSF                     | 6.06  | 10.30                        | 5.64                                      | 8.41                         |
| sCD137/4-1BB/<br>TNFRSF9  | 5.42  | 10.47                        | 6.42                                      | 15.08                        |
| sICAM-1                   | 78.00   | 135.14                       | 33.35                                     | 63.28                        |
| CXCL16                    | 5.91  | 11.10                        | 5.82                                      | 13.15                        |
| MCP-5                     | 0.68  | 1.10                         | 0.81                                      | 1.35                         |
| MIP-2                     | 0.26  | 0.57                         | 0.29                                      | 0.72                         |
| MIG                       | 15.57   | 24.26                        | 13.97                                     | 21.56                        |
| RANTES                    | 6.10  | 9.22                         | 5.29                                      | 11.76                        |
| VEGF-A                    | 3.12  | 4.06                         | 2.08                                      | 3.56                         |
| TNF $\alpha$              | 2.07  | 3.72                         | 1.71                                      | 4.29                         |
| IL-11                     | 13.06   | 15.87                        | 9.51                                      | 10.25                        |

For research use only. Not for use in diagnostic procedures.

| <b>Analyte</b>     | <b>Overnight Protocol<br/>(n = 13 Assays)</b> |                              | <b>2-Hour Protocol<br/>(n = 3 Assays)</b> |                              |
|--------------------|---|------------------------------|---|------------------------------|
|                    | <b>MinDC<br/>(pg/mL)</b>                      | <b>MinDC+2SD<br/>(pg/mL)</b> | <b>MinDC<br/>(pg/mL)</b>                  | <b>MinDC+2SD<br/>(pg/mL)</b> |
| sRAGE              | 24.5  | 49.43                        | 21.82                                     | 44.58                        |
| CCL25/TECK         | 15.51   | 27.79                        | 18.28                                     | 30.06                        |
| BAFF               | 1.76  | 3.66                         | 3.52                                      | 10.53                        |
| GDF-15             | 0.55  | 0.98                         | 0.81                                      | 1.84                         |
| FLT3L              | 2.57  | 4.43                         | 1.23                                      | 1.73                         |
| CHI3L1             | 17.23   | 24.34                        | 7.84                                      | 13.30                        |
| MCP-2              | 0.27  | 0.51                         | 0.32                                      | 0.94                         |
| MCP-3              | 1.28  | 1.71                         | 1.38                                      | 1.75                         |
| Erythropoietin     | 8.39  | 13.77                        | 6.07                                      | 15.55                        |
| Fractalkine/CX3CL1 | 0.67  | 1.24                         | 0.49                                      | 0.64                         |
| IL-31              | 32.68   | 60.12                        | 32.83                                     | 76.18                        |
| Betacellulin       | 1.46  | 2.86                         | 1.95                                      | 4.34                         |

## Precision

Intra-assay precision is generated from the mean of the % CVs from 8 reportable results across two different concentrations of analytes in a single assay. Inter-assay precision is generated from the mean of the % CV's across two different concentrations of analytes across 8 different assays for overnight and 3 different assays for same day (see next page).

| Analyte                | Overnight Protocol |                 | Same Day Protocol |                 |
|------------------------|--------------------|-----------------|-------------------|-----------------|
|                        | Inter-assay %CV    | Intra-assay %CV | Inter-assay %CV   | Intra-assay %CV |
| sFasL                  | <15%               | <10%            | <15%              | <10%            |
| IL-22                  | <15%               | <10%            | <15%              | <10%            |
| G-CSF                  | <15%               | <10%            | <15%              | <10%            |
| Eotaxin/CCL11          | <15%               | <10%            | <15%              | <10%            |
| GM-CSF                 | <15%               | <10%            | <15%              | <10%            |
| IL-33                  | <15%               | <10%            | <15%              | <10%            |
| IFN $\gamma$           | <15%               | <10%            | <15%              | <10%            |
| IL-1 $\alpha$          | <15%               | <10%            | <15%              | <10%            |
| IL-17F                 | <15%               | <10%            | <15%              | <10%            |
| IL-1 $\beta$           | <15%               | <10%            | <15%              | <10%            |
| IL-2                   | <15%               | <10%            | <15%              | <10%            |
| IFN $\beta$            | <15%               | <10%            | <15%              | <10%            |
| IL-4                   | <15%               | <10%            | <15%              | <10%            |
| IL-3                   | <15%               | <10%            | <15%              | <10%            |
| IL-5                   | <15%               | <10%            | <15%              | <10%            |
| IL-18                  | <15%               | <10%            | <15%              | <10%            |
| IL-6                   | <15%               | <10%            | <15%              | <10%            |
| FGF-2                  | <15%               | <10%            | <15%              | <10%            |
| IL-7                   | <15%               | <10%            | <15%              | <10%            |
| EGF                    | <15%               | <10%            | <15%              | <10%            |
| IL-9                   | <15%               | <10%            | <15%              | <10%            |
| IFN $\alpha$           | <15%               | <10%            | <15%              | <10%            |
| IL-28B/IFN $\lambda$ 3 | <15%               | <10%            | <15%              | <10%            |
| IL-10                  | <15%               | <10%            | <15%              | <10%            |
| MIP-3 $\alpha$         | <15%               | <10%            | <15%              | <10%            |
| IL-12p40               | <15%               | <10%            | <15%              | <10%            |
| Granzyme B             | <15%               | <10%            | <15%              | <10%            |
| IL-12p70               | <15%               | <10%            | <15%              | <10%            |

For research use only. Not for use in diagnostic procedures.

| <b>Analyte</b>            | <b>Overnight Protocol</b> |                        | <b>Same Day Protocol</b> |                        |
|---------------------------|---------------------------|------------------------|--------------------------|------------------------|
|                           | <b>Inter-assay %CV</b>    | <b>Intra-assay %CV</b> | <b>Inter-assay %CV</b>   | <b>Intra-assay %CV</b> |
| TARC                      | <15%                      | <10%                   | <15%                     | <10%                   |
| IL-16                     | <15%                      | <10%                   | <15%                     | <10%                   |
| LIF                       | <15%                      | <10%                   | <15%                     | <10%                   |
| IL-13                     | <15%                      | <10%                   | <15%                     | <10%                   |
| LIX                       | <15%                      | <10%                   | <15%                     | <10%                   |
| IL-15                     | <15%                      | <10%                   | <15%                     | <10%                   |
| Exodus-2/<br>CCL21/6Ckine | <20%                      | <10%                   | <20%                     | <10%                   |
| IL-17A                    | <15%                      | <10%                   | <15%                     | <10%                   |
| IP-10                     | <15%                      | <10%                   | <15%                     | <10%                   |
| Fas                       | <15%                      | <10%                   | <15%                     | <10%                   |
| IL-20                     | <15%                      | <10%                   | <15%                     | <10%                   |
| KC                        | <15%                      | <10%                   | <15%                     | <10%                   |
| MCP-1                     | <15%                      | <10%                   | <15%                     | <10%                   |
| MIP-3 $\beta$             | <15%                      | <10%                   | <15%                     | <10%                   |
| MIP-1 $\alpha$            | <15%                      | <10%                   | <15%                     | <10%                   |
| MDC                       | <15%                      | <10%                   | <15%                     | <10%                   |
| MIP-1 $\beta$             | <15%                      | <10%                   | <15%                     | <10%                   |
| M-CSF                     | <15%                      | <10%                   | <15%                     | <10%                   |
| sCD137/4-1BB/<br>TNFRSF9  | <15%                      | <10%                   | <15%                     | <10%                   |
| sICAM-1                   | <15%                      | <10%                   | <15%                     | <10%                   |
| CXCL16                    | <15%                      | <10%                   | <15%                     | <10%                   |
| MCP-5                     | <15%                      | <10%                   | <15%                     | <10%                   |
| MIP-2                     | <15%                      | <10%                   | <15%                     | <10%                   |
| MIG                       | <15%                      | <10%                   | <15%                     | <10%                   |
| RANTES                    | <15%                      | <10%                   | <15%                     | <10%                   |
| VEGF-A                    | <15%                      | <10%                   | <15%                     | <10%                   |
| TNF $\alpha$              | <15%                      | <10%                   | <15%                     | <10%                   |

For research use only. Not for use in diagnostic procedures.

| Analyte            | Overnight Protocol |                 | Same Day Protocol |                 |
|--------------------|--------------------|-----------------|-------------------|-----------------|
|                    | Inter-assay %CV    | Intra-assay %CV | Inter-assay %CV   | Intra-assay %CV |
| IL-11              | <15%               | <10%            | <15%              | <10%            |
| sRAGE              | <15%               | <10%            | <15%              | <10%            |
| CCL25/TECK         | <15%               | <10%            | <15%              | <10%            |
| BAFF               | <15%               | <10%            | <15%              | <10%            |
| GDF-15             | <15%               | <10%            | <15%              | <10%            |
| FLT3L              | <15%               | <10%            | <15%              | <10%            |
| CHI3L1             | <15%               | <10%            | <15%              | <10%            |
| MCP-2              | <15%               | <10%            | <15%              | <10%            |
| MCP-3              | <15%               | <10%            | <15%              | <10%            |
| Erythropoietin     | <15%               | <10%            | <15%              | <10%            |
| Fractalkine/CX3CL1 | <15%               | <10%            | <15%              | <10%            |
| IL-31              | <15%               | <10%            | <15%              | <10%            |
| Betacellulin       | <15%               | <10%            | <15%              | <10%            |

## Accuracy

**Spike Recovery:** The data represent mean percent recovery of spiked standards ranging from low, medium, and high concentration in serum matrices (n = 5, see next page).

| <b>Analyte</b>         | <b>Overnight Protocol</b>         | <b>2-Hour Protocol</b>            |
|------------------------|-----------------------------------|-----------------------------------|
|                        | <b>% Recovery in Serum Matrix</b> | <b>% Recovery in Serum Matrix</b> |
| sFasL                  | 94                                | 103                               |
| IL-22                  | 90                                | 96                                |
| G-CSF                  | 90                                | 110                               |
| Eotaxin/CCL11          | 88                                | 100                               |
| GM-CSF                 | 98                                | 97                                |
| IL-33                  | 93                                | 99                                |
| IFN $\gamma$           | 94                                | 100                               |
| IL-1 $\alpha$          | 94                                | 94                                |
| IL-17F                 | 91                                | 103                               |
| IL-1 $\beta$           | 93                                | 102                               |
| IL-2                   | 83                                | 89                                |
| IFN $\beta$            | 95                                | 100                               |
| IL-4                   | 89                                | 92                                |
| IL-3                   | 94                                | 98                                |
| IL-5                   | 83                                | 95                                |
| IL-18                  | 87                                | 84                                |
| IL-6                   | 89                                | 97                                |
| FGF-2                  | 96                                | 96                                |
| IL-7                   | 94                                | 98                                |
| EGF                    | 97                                | 104                               |
| IL-9                   | 93                                | 96                                |
| IFN $\alpha$           | 91                                | 98                                |
| IL-28B/IFN $\lambda$ 3 | 89                                | 105                               |
| IL-10                  | 93                                | 100                               |
| MIP-3 $\alpha$         | 95                                | 103                               |
| IL-12p40               | 91                                | 96                                |
| Granzyme B             | 93                                | 109                               |
| IL-12p70               | 92                                | 91                                |

For research use only. Not for use in diagnostic procedures.



| <b>Analyte</b>        | <b>Overnight Protocol</b>         | <b>2-Hour Protocol</b>            |
|-----------------------|-----------------------------------|-----------------------------------|
|                       | <b>% Recovery in Serum Matrix</b> | <b>% Recovery in Serum Matrix</b> |
| TARC                  | 91                                | 102                               |
| IL-16                 | 95                                | 101                               |
| LIF                   | 90                                | 101                               |
| IL-13                 | 92                                | 99                                |
| LIX                   | 92                                | 99                                |
| IL-15                 | 88                                | 102                               |
| Exodus-2/CCL21/6Ckine | 100                               | 110                               |
| IL-17A                | 89                                | 92                                |
| IP-10                 | 92                                | 103                               |
| Fas                   | 92                                | 102                               |
| IL-20                 | 99                                | 94                                |
| KC                    | 88                                | 103                               |
| MCP-1                 | 92                                | 93                                |
| MIP-3 $\beta$         | 93                                | 101                               |
| MIP-1 $\alpha$        | 97                                | 93                                |
| MDC                   | 90                                | 96                                |
| MIP-1 $\beta$         | 97                                | 94                                |
| M-CSF                 | 93                                | 98                                |
| sCD137/4-1BB/TNFRSF9  | 95                                | 104                               |
| sICAM-1               | 89                                | 90                                |
| CXCL16                | 94                                | 106                               |
| MCP-5                 | 91                                | 103                               |
| MIP-2                 | 87                                | 92                                |
| MIG                   | 93                                | 99                                |
| RANTES                | 93                                | 97                                |
| VEGF-A                | 94                                | 108                               |
| TNF $\alpha$          | 90                                | 92                                |
| IL-11                 | 90                                | 94                                |
| sRAGE                 | 92                                | 99                                |

For research use only. Not for use in diagnostic procedures.

| <b>Analyte</b>     | <b>Overnight Protocol</b>         | <b>2-Hour Protocol</b>            |
|--------------------|-----------------------------------|-----------------------------------|
|                    | <b>% Recovery in Serum Matrix</b> | <b>% Recovery in Serum Matrix</b> |
| CCL25/TECK         | 91                                | 104                               |
| BAFF               | 91                                | 102                               |
| GDF-15             | 95                                | 99                                |
| FLT3L              | 94                                | 99                                |
| CHI3L1             | 95                                | 102                               |
| MCP-2              | 92                                | 97                                |
| MCP-3              | 91                                | 100                               |
| Erythropoietin     | 93                                | 102                               |
| Fractalkine/CX3CL1 | 92                                | 104                               |
| IL-31              | 92                                | 100                               |
| Betacellulin       | 96                                | 106                               |

## Troubleshooting

| <b>Problem</b>          | <b>Probable Cause</b>   | <b>Solution</b>  |
|-------------------------|---|--|
|                         | Plate washer aspirate height set too low                          | Adjust aspiration height according to manufacturers' instructions.   |
|                         | Bead mix prepared inappropriately                                 | Sonicate bead vials and vortex just prior to adding to bead mix bottle according to protocol. Agitate bead mix intermittently in reservoir while pipetting this into the plate.  |
|                         | Samples cause interference due to particulate matter or viscosity | See above. Also sample probe may need to be cleaned with alcohol flushes, back flushes, and washes; or, if needed, probe should be removed and sonicated.  |
| Insufficient bead count | Probe height not adjusted correctly                               | <p>When reading the assay on Luminex® 200™, adjust probe height to the kit solid plate or to the recommended filter plates using 3 alignment discs. When reading the assay on MAGPIX®, adjust probe height to the kit solid plate or to the recommended filter plates using 2 alignment discs. When reading the assay on FLEXMAP 3D®, adjust probe height to the kit solid plate using 1 alignment disc. For FLEXMAP 3D® when using the solid plate in the kit, the final resuspension should be with 150 µL Sheath Fluid PLUS in each well and 75 µL should be aspirated.</p> <p>When reading the assay on xMAP® INTELLIFLEX, adjust probe height based on the type of plate you are using, place an alignment disk or an alignment sphere in the well according to the protocol recommended by Luminex®.</p> |

| <b>Problem</b>              | <b>Probable Cause</b>                                    | <b>Solution</b>  |
|-----------------------------|--|--|
| Background is too high      | Background wells were contaminated                       | Avoid cross-well contamination by using sealer appropriately and pipetting with multichannel pipettes without touching reagent in plate.                           |
|                             | Matrix used has endogenous analyte or interference       | Check matrix ingredients for cross-reacting components (for example, interleukin modified tissue culture medium).  |
|                             | Insufficient washes                                      | Increase number of washes.   |
| Beads not in region or gate | Luminex® instrument not calibrated correctly or recently | Calibrate Luminex® instrument based on manufacturer's instructions, at least once a week or if temperature has changed by >3 °C.                                   |
|                             | Gate settings not adjusted correctly                     | Some Luminex® instruments (for example Bio-Plex®) require different gate settings than those described in the kit protocol. Use instrument default settings.       |
|                             | Wrong bead regions in protocol template                  | Check kit protocol for correct bead regions or analyte selection.  |
|                             | Incorrect sample type used                               | Samples containing organic solvents or if highly viscous should be diluted or dialyzed as required.  |
|                             | Instrument not washed or primed                          | Prime the Luminex® instrument 4 times to rid it of air bubbles, wash 4 times with Sheath Fluid PLUS or water if there is any remnant alcohol or sanitizing liquid. |
|                             | Beads were exposed to light                              | Keep plate and bead mix covered with dark lid or aluminum foil during all incubation steps.  |

| <b>Problem</b>                                  | <b>Probable Cause</b>   | <b>Solution</b>   |
|---|---|---|
| Signal for whole plate is same as background    | Incorrect or no Detection Antibody was added  | Add appropriate Detection Antibody and continue.  |
|   | Streptavidin-Phycoerythrin was not added  | Add Streptavidin-Phycoerythrin according to protocol. If Detection Antibody has already been removed, sensitivity may be low.   |
| Low signal for standard curve                   | Detection Antibody may have been removed prior to adding Streptavidin-Phycoerythrin | May need to repeat assay if desired sensitivity not achieved.   |
|   | Incubations done at inappropriate temperatures, timings, or agitation               | Assay conditions need to be checked.  |
| Signals too high, standard curves are saturated | Calibration target value set too high   | With some Luminex® instruments (for example, Bio-Plex®) default target setting for RP1 calibrator is set at high PMT. Use low target value for calibration and reanalyze plate. |
|   | Plate incubation was too long with standard curve and samples                       | Use shorter incubation time.  |
| Sample readings are out of range                | Samples contain no or below detectable levels of analyte                            | If below detectable levels, it may be possible to use higher sample volume. Check with technical support for appropriate protocol modifications.                                |
|   | Samples contain analyte concentrations higher than highest standard point           | Samples may require dilution and reanalysis for just that particular analyte.   |
|   | Standard curve was saturated at higher end of curve                                 | See above.  |

For research use only. Not for use in diagnostic procedures.

| <b>Problem</b>                             | <b>Probable Cause</b>  | <b>Solution</b>   |
|--|--|---|
| High variation in samples and/or standards | Multichannel pipette may not be calibrated                               | Calibrate pipettes.   |
|  | Plate washing was not uniform  | Confirm all reagents are removed completely in all wash steps.  |
|  | Samples may have high particulate matter or other interfering substances | See above.  |
|  | Plate agitation was insufficient   | Plate should be agitated during all incubation steps using an orbital plate shaker at a speed where beads are in constant motion without causing splashing.   |
|  | Cross-well contamination   | Check when reusing plate sealer that no reagent has touched sealer. Care should be taken when using same pipette tips that are used for reagent additions and that pipette tip does not touch reagent in plate. |

## Product Ordering

Order products online at [SigmaAldrich.com](http://SigmaAldrich.com).

| Replacement Reagents   | Note                                      | Catalogue No.    |
|--|---|------------------|
| Mouse Cytokine Expansion Panel 1 Standard 1                  | For configurable kit                      | MCYT1-8190-1     |
| Mouse Cytokine Expansion Panel 1 Standard 2                  | For 41-plex, 68-plex, or configurable kit | MCYT1-8190-2     |
| Mouse Cytokine Expansion Panel 1 QC 1 & 2 for Standard 1     | For configurable kit                      | MCYT1-6190-1     |
| Mouse Cytokine Expansion Panel 1 QC 1 & 2 for Standard 2     | For 41-plex, 68-plex, or configurable kit | MCYT1-6190-2     |
| Serum Matrix   | -   | MXMSM-MCYT1      |
| Mouse Cytokine Expansion Panel 1 Detection Antibodies 1      | For configurable kit                      | MCYT1-1190-1     |
| Mouse Cytokine Expansion Panel 1 Detection Antibodies 2      | For 41-plex, 68-plex, or configurable kit | MCYT1-1190-2     |
| Streptavidin-Phycoerythrin                                   | -   | L-SAPE20         |
| Assay Buffer   | -   | L-AB             |
| Set of two 96-Well plates with sealers                       | -   | MAG-PLATE        |
| 10X Wash Buffer  | -   | L-WB             |
| Mouse Cytokine Expansion Panel 1 S-Plex Premixed Beads*      | -   | MCYT1PMXS-MG     |
| Mouse Cytokine Expansion Panel 1 L-Plex Premixed Beads*      | -   | MCYT1PMXL-MG     |
| MILLIPLEX® Mouse Cytokine/Chemokine/GF Exp Pnl 1 Sm-Plex PMX | -   | MCYT1-190K-SPX   |
| MILLIPLEX® Mouse Cytokine/ChemokineGF Exp Pnl 1 L-Plex PMX   | -   | MCYT1-190K-LPX   |
| MILLIPLEX® Mu Cytokine/Chemokine/GF Exp Pnl 1 BK Sm-Plex PMX | BULK PACKAGING                            | MCYT1-190K-SPXBK |
| MILLIPLEX® Mu Cytokine/Chemokine/GF Exp Pnl 1 BLK L-Plex PMX | BULK PACKAGING                            | MCYT1-190K-LPXBK |

\* For individual beads, see next page.

For research use only. Not for use in diagnostic procedures.

## Antibody-Immobilized Magnetic Beads

| <b>Bead/Analyte Name</b>               | <b>Bead No.</b> | <b>Catalogue No</b> |
|--|-----------------|---------------------|
| Anti-Mouse sFasL Bead                  | 9               | MSFASL-MG           |
| Anti-Mouse IL-22 Bead                  | 12              | MIL22-MG            |
| Anti-Mouse G-CSF Bead                  | 13              | MGCSF-MG            |
| Anti-Mouse Eotaxin/CCL11 Bead          | 14              | MCCL11-MG           |
| Anti-Mouse GM-CSF Bead                 | 15              | MGMCSF-MG           |
| Anti-Mouse IL-33 Bead                  | 18              | MIL33-MG            |
| Anti-Mouse IFN $\gamma$ Bead           | 19              | MIFNY-MG            |
| Anti-Mouse IL-1 $\alpha$ Bead          | 21              | MIL1A-MG            |
| Anti-Mouse IL-17F Bead                 | 22              | MIL17F-MG           |
| Anti-Mouse IL-1 $\beta$ Bead           | 25              | MIL1B-MG            |
| Anti- Mouse IL-2 Bead                  | 26              | MIL2-MG             |
| Anti-Mouse IFN $\beta$ Bead            | 27              | MIFNB-MG            |
| Anti-Mouse IL-4 Bead                   | 28              | MIL4-MG             |
| Anti-Mouse IL-3 Bead                   | 29              | MIL3-MG             |
| Anti-Mouse IL-5 Bead                   | 30              | MIL5-MG             |
| Anti-Mouse IL-18 Bead                  | 33              | MIL18-MG            |
| Anti-Mouse IL-6 Bead                   | 34              | MIL6-MG             |
| Anti-Mouse FGF-2 Bead                  | 35              | MFGF2-MG            |
| Anti-Mouse IL-7 Bead                   | 36              | MIL7-MG             |
| Anti-Mouse EGF Bead                    | 37              | MEGF-MG             |
| Anti-Mouse IL-9 Bead                   | 38              | MIL9-MG             |
| Anti-Mouse IFN $\alpha$ Bead           | 39              | MIFNA-MG            |
| Anti-Mouse IL-28B/IFN $\lambda$ 3 Bead | 42              | MIL28B-MG           |
| Anti-Mouse IL-10 Bead                  | 43              | MIL10-MG            |
| Anti-Mouse MIP-3 $\alpha$ Bead         | 44              | MMIP3A-MG           |
| Anti-Mouse IL-12p40 Bead               | 45              | MIL12P40-MG         |
| Anti-Mouse Granzyme B Bead             | 46              | MGZMB-MG            |

For research use only. Not for use in diagnostic procedures.



| <b>Bead/Analyte Name</b>                  | <b>Bead No.</b> | <b>Catalogue No</b> |
|---|-----------------|---------------------|
| Anti-Mouse IL-12p70 Bead                  | 47              | MIL12P70-MG         |
| Anti-Mouse TARC Bead                      | 48              | MTARC-MG            |
| Anti-Mouse IL-16 Bead                     | 49              | MIL16-MG            |
| Anti-Mouse LIF Bead                       | 51              | MLIF-MG             |
| Anti-Mouse IL-13 Bead                     | 52              | MIL13-MG            |
| Anti-Mouse LIX Bead                       | 53              | MLIX-MG             |
| Anti-Mouse IL-15 Bead                     | 54              | MIL15-MG            |
| Anti-Mouse Exodus-2/<br>CCL21/6Ckine Bead | 55              | MCCL21-MG           |
| Anti-Mouse IL-17A Bead                    | 56              | MIL17A-MG           |
| Anti-Mouse IP-10 Bead                     | 57              | MIP10-MG            |
| Anti-Mouse Fas Bead                       | 58              | MFAS-MG             |
| Anti-Mouse IL-20 Bead                     | 59              | MIL20-MG            |
| Anti-Mouse KC Bead                        | 61              | MKC-MG              |
| Anti-Mouse MCP-1 Bead                     | 62              | MMCP1-MG            |
| Anti-Mouse MIP-3 $\beta$ Bead             | 63              | MMIP3B-MG           |
| Anti-Mouse MIP-1 $\alpha$ Bead            | 64              | MMIP1A-MG           |
| Anti-Mouse MDC Bead                       | 65              | MMDC-MG             |
| Anti-Mouse MIP-1 beta Bead                | 66              | MMIP1B-MG           |
| Anti-Mouse M-CSF Bead                     | 67              | MMCSF-MG            |
| Anti-Mouse sCD137/4-1BB/<br>TNFRSF9 Bead  | 68              | MCD137-MG           |
| Anti-Mouse sICAM-1 Bead                   | 69              | MSICAM1-MG          |
| Anti-Mouse CXCL16 Bead                    | 70              | MCXCL16-MG          |
| Anti-Mouse MCP-5 Bead                     | 72              | MMCP5-MG            |
| Anti-Mouse MIP-2 Bead                     | 73              | MMIP2-MG            |
| Anti-Mouse MIG Bead                       | 74              | MMIG-MG             |
| Anti-Mouse RANTES Bead                    | 75              | MRANTES-MG          |
| Anti-Mouse VEGF-A Bead                    | 76              | MVEGFA-MG           |

For research use only. Not for use in diagnostic procedures.

| <b>Bead/Analyte Name</b>               | <b>Bead No.</b> | <b>Catalogue No</b> |
|--|-----------------|---------------------|
| Anti-Mouse TNF $\alpha$ Bead           | 77              | MTNFA-MG            |
| Anti-Mouse IL-11 Bead                  | 78              | MIL11-MG            |
| Anti-Mouse sRAGE Bead                  | 79              | MSRAGE-MG           |
| Anti-Mouse CCL25/TECK Bead             | 80              | MCCL25-MG           |
| Anti-Mouse BAFF Bead                   | 81              | MBAFF-MG            |
| Anti-Mouse GDF-15 Bead                 | 82              | MGDF15-MG           |
| Anti-Mouse FLT3L Bead                  | 83              | MFLT3L-MG           |
| Anti-Mouse CHI3L1 Bead                 | 84              | MCHI3L1-MG          |
| Anti-Mouse MCP-2 Bead                  | 85              | MMCP2-MG            |
| Anti-Mouse MCP-3 Bead                  | 86              | MMCP3-MG            |
| Anti-Mouse Erythropoietin Bead         | 87              | MEPO-MG             |
| Anti-Mouse Fractalkine/<br>CX3CL1 Bead | 88              | MCX3CL1-MG          |
| Anti-Mouse IL-31 Bead                  | 89              | MIL31-MG            |
| Anti-Mouse Betacellulin Bead           | 90              | MBTC-MG             |

| Analyte/<br>Bead Name | Luminex® Magnetic<br>Bead Region | Customizable 68 Analytes<br>(70X concentration, 70 µL) |                | MCYT1PMXS-MG<br>(41-Plex Premixed Beads) | MCYT1PMXL-MG<br>(68-Plex Premixed Beads) | MCYT1-8190-1<br>(35-plex Standard Mix) | MCYT1-8190-2<br>(68-plex Standard Mix) | MCYT1-1190-1<br>(35-plex Detection Mix) | MCYT1-1190-2<br>(68-plex Detection Mix) |
|-----------------------|----------------------------------|--|----------------|--|--|--|--|---|---|
|                       |                                  | Available  | Catalog Number |  |  |  |  |   |   |
| sFasL                 | 9                                | ✓  | MSFASL-MG      |  | ✓  |  | ✓                                      |   | ✓                                       |
| IL-22                 | 12                               | ✓  | MIL22-MG       | ✓  | ✓  |  | ✓                                      |   | ✓                                       |
| G-CSF                 | 13                               | ✓  | MGCSF-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| Eotaxin/CCL11         | 14                               | ✓  | MCCL11-MG      | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| GM-CSF                | 15                               | ✓  | MGMCSF-MG      | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IL-33                 | 18                               | ✓  | MIL33-MG       | ✓  | ✓  |  | ✓                                      |   | ✓                                       |
| IFN $\gamma$          | 19                               | ✓  | MIFNY-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IL-1 $\alpha$         | 21                               | ✓  | MIL1A-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IL-17F                | 22                               | ✓  | MIL17F-MG      | ✓  | ✓  |  | ✓                                      |   | ✓                                       |
| IL-1 $\beta$          | 25                               | ✓  | MIL1B-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IL-2                  | 26                               | ✓  | MIL2-MG        | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IFN $\beta$           | 27                               | ✓  | MIFNB-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IL-4                  | 28                               | ✓  | MIL4-MG        | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IL-3                  | 29                               | ✓  | MIL3-MG        | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IL-5                  | 30                               | ✓  | MIL5-MG        | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IL-18                 | 33                               | ✓  | MIL18-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IL-6                  | 34                               | ✓  | MIL6-MG        | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| FGF-2                 | 35                               | ✓  | MFGF2-MG       | ✓  | ✓  |  | ✓                                      |   | ✓                                       |

For research use only. Not for use in diagnostic procedures.

| Analyte/<br>Bead Name     | Luminex® Magnetic<br>Bead Region | Customizable 68 Analytes<br>(70X concentration, 70 µL) |                | MCYT1PMXS-MG<br>(41-Plex Premixed Beads) | MCYT1PMXL-MG<br>(68-Plex Premixed Beads) | MCYT1-8190-1<br>(35-plex Standard Mix) | MCYT1-8190-2<br>(68-plex Standard Mix) | MCYT1-1190-1<br>(35-plex Detection Mix) | MCYT1-1190-2<br>(68-plex Detection Mix) |
|---------------------------|----------------------------------|--|----------------|--|--|--|--|---|---|
|                           |                                  | Available  | Catalog Number |  |  |  |  |   |   |
| IL-7                      | 36                               | ✓  | MIL7-MG        | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| EGF                       | 37                               | ✓  | MEGF-MG        | ✓  | ✓  |  | ✓                                      |   | ✓                                       |
| IL-9                      | 38                               | ✓  | MIL9-MG        | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IFNα                      | 39                               | ✓  | MIFNA-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IL-28B/IFNλ3              | 42                               | ✓  | MIL28B-MG      | ✓  | ✓  |  | ✓                                      |   | ✓                                       |
| IL-10                     | 43                               | ✓  | MIL10-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| MIP-3α                    | 44                               | ✓  | MMIP3A-MG      |  | ✓  |  | ✓                                      |   | ✓                                       |
| IL-12p40                  | 45                               | ✓  | MIL12P40-MG    | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| Granzyme B                | 46                               | ✓  | MGZMB-MG       |  | ✓  |  | ✓                                      |   | ✓                                       |
| IL-12p70                  | 47                               | ✓  | MIL12P70-MG    | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| TARC                      | 48                               | ✓  | MTARC-MG       |  | ✓  |  | ✓                                      |   | ✓                                       |
| IL-16                     | 49                               | ✓  | MIL16-MG       |  | ✓  |  | ✓                                      |   | ✓                                       |
| LIF                       | 51                               | ✓  | MLIF-MG        | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IL-13                     | 52                               | ✓  | MIL13-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| LIX                       | 53                               | ✓  | MLIX-MG        | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IL-15                     | 54                               | ✓  | MIL15-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| Exodus-2/<br>CCL21/6Ckine | 55                               | ✓  | MCCL21-MG      |  | ✓  |  | ✓                                      |   | ✓                                       |
| IL-17A                    | 56                               | ✓  | MIL17A-MG      | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |

For research use only. Not for use in diagnostic procedures.

| Analyte/<br>Bead Name    | Luminex® Magnetic<br>Bead Region | Customizable 68 Analytes<br>(70X concentration, 70 µL) |                | MCYT1PMXS-MG<br>(41-Plex Premixed Beads) | MCYT1PMXL-MG<br>(68-Plex Premixed Beads) | MCYT1-8190-1<br>(35-plex Standard Mix) | MCYT1-8190-2<br>(68-plex Standard Mix) | MCYT1-1190-1<br>(35-plex Detection Mix) | MCYT1-1190-2<br>(68-plex Detection Mix) |
|--------------------------|----------------------------------|--|----------------|--|--|--|--|---|---|
|                          |                                  | Available  | Catalog Number |  |  |  |  |   |   |
| IP-10                    | 57                               | ✓  | MIP10-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| Fas                      | 58                               | ✓  | MFAS-MG        |  | ✓  |  | ✓                                      |   | ✓                                       |
| IL-20                    | 59                               | ✓  | MIL20-MG       |  | ✓  |  | ✓                                      |   | ✓                                       |
| KC                       | 61                               | ✓  | MKC-MG         | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| MCP-1                    | 62                               | ✓  | MMCP1-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| MIP-3β                   | 63                               | ✓  | MMIP3B-MG      |  | ✓  |  | ✓                                      |   | ✓                                       |
| MIP-1α                   | 64                               | ✓  | MMIP1A-MG      | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| MDC                      | 65                               | ✓  | MMDC-MG        |  | ✓  |  | ✓                                      |   | ✓                                       |
| MIP-1β                   | 66                               | ✓  | MMIP1B-MG      | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| M-CSF                    | 67                               | ✓  | MMCSF-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| sCD137/4-1BB/<br>TNFRSF9 | 68                               | ✓  | MCD137-MG      |  | ✓  |  | ✓                                      |   | ✓                                       |
| sICAM-1                  | 69                               | ✓  | MSICAM1-MG     |  | ✓  |  | ✓                                      |   | ✓                                       |
| CXCL16                   | 70                               | ✓  | MCXCL16-MG     |  | ✓  |  | ✓                                      |   | ✓                                       |
| MCP-5                    | 72                               | ✓  | MMCP5-MG       |  | ✓  |  | ✓                                      |   | ✓                                       |
| MIP-2                    | 73                               | ✓  | MMIP2-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| MIG                      | 74                               | ✓  | MMIG-MG        | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| RANTES                   | 75                               | ✓  | MRANTES-MG     | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| VEGF-A                   | 76                               | ✓  | MVEGFA-MG      | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |

For research use only. Not for use in diagnostic procedures.

| Analyte/<br>Bead Name  | Luminex® Magnetic<br>Bead Region | Customizable 68 Analytes<br>(70X concentration, 70 µL) |                | MCYT1PMXS-MG<br>(41-Plex Premixed Beads) | MCYT1PMXL-MG<br>(68-Plex Premixed Beads) | MCYT1-8190-1<br>(35-plex Standard Mix) | MCYT1-8190-2<br>(68-plex Standard Mix) | MCYT1-1190-1<br>(35-plex Detection Mix) | MCYT1-1190-2<br>(68-plex Detection Mix) |
|------------------------|----------------------------------|--|----------------|--|--|--|--|---|---|
|                        |                                  | Available  | Catalog Number |  |  |  |  |   |   |
| TNFα                   | 77                               | ✓  | MTNFA-MG       | ✓  | ✓  | ✓                                      | ✓                                      | ✓                                       | ✓                                       |
| IL-11                  | 78                               | ✓  | MIL11-MG       |  | ✓  |  | ✓                                      |   | ✓                                       |
| sRAGE                  | 79                               | ✓  | MSRAGE-MG      |  | ✓  |  | ✓                                      |   | ✓                                       |
| CCL25/TECK             | 80                               | ✓  | MCCL25-MG      |  | ✓  |  | ✓                                      |   | ✓                                       |
| BAFF                   | 81                               | ✓  | MBAFF-MG       |  | ✓  |  | ✓                                      |   | ✓                                       |
| GDF-15                 | 82                               | ✓  | MGDF15-MG      |  | ✓  |  | ✓                                      |   | ✓                                       |
| FLT3L                  | 83                               | ✓  | MFLT3L-MG      |  | ✓  |  | ✓                                      |   | ✓                                       |
| CHI3L1                 | 84                               | ✓  | MCHI3L1-MG     |  | ✓  |  | ✓                                      |   | ✓                                       |
| MCP-2                  | 85                               | ✓  | MMCP2-MG       |  | ✓  |  | ✓                                      |   | ✓                                       |
| MCP-3                  | 86                               | ✓  | MMCP3-MG       |  | ✓  |  | ✓                                      |   | ✓                                       |
| Erythropoietin         | 87                               | ✓  | MEPO-MG        |  | ✓  |  | ✓                                      |   | ✓                                       |
| Fractalkine/<br>CX3CL1 | 88                               | ✓  | MCX3CL1-MG     |  | ✓  |  | ✓                                      |   | ✓                                       |
| IL-31                  | 89                               | ✓  | MIL31-MG       |  | ✓  |  | ✓                                      |   | ✓                                       |
| Betacellulin           | 90                               | ✓  | MBTC-MG        |  | ✓  |  | ✓                                      |   | ✓                                       |

For research use only. Not for use in diagnostic procedures.

## Well Map

|   | 1                             | 2          | 3               | 4    | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
|---|-------------------------------|------------|-----------------|------|---|---|---|---|---|----|----|----|
| A | 0<br>Standard<br>(Background) | Standard 4 | QC-1<br>Control | Etc. |   |   |   |   |   |    |    |    |
| B | 0<br>Standard<br>(Background) | Standard 4 | QC-1<br>Control |      |   |   |   |   |   |    |    |    |
| C | Standard 1                    | Standard 5 | QC-2<br>Control |      |   |   |   |   |   |    |    |    |
| D | Standard 1                    | Standard 5 | QC-2<br>Control |      |   |   |   |   |   |    |    |    |
| E | Standard 2                    | Standard 6 | Sample<br>1     |      |   |   |   |   |   |    |    |    |
| F | Standard 2                    | Standard 6 | Sample<br>1     |      |   |   |   |   |   |    |    |    |
| G | Standard 3                    | Standard 7 | Sample<br>2     |      |   |   |   |   |   |    |    |    |
| H | Standard 3                    | Standard 7 | Sample<br>2     |      |   |   |   |   |   |    |    |    |

---

## Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

### Technical Assistance

Visit the tech service page at [SigmaAldrich.com/techservice](https://SigmaAldrich.com/techservice).

### Standard Warranty

The applicable warranty for the products listed in this publication may be found at [SigmaAldrich.com/terms](https://SigmaAldrich.com/terms).

### Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://SigmaAldrich.com/offices).

### Safety Data Sheets (SDS)

Safety Data Sheets are available on the product page at [SigmaAldrich.com](https://SigmaAldrich.com).

The life science business of Merck KGaA, Darmstadt, Germany operates as MilliporeSigma in the U.S. and Canada.

MilliporeSigma, Millipore, Milliplex and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

© 2025 Merck KGaA, Darmstadt, Germany and/or its affiliates.  
All Rights Reserved.

For research use only. Not for use in diagnostic procedures.

23166000 Rev 01/25

The logo for MilliporeSigma, featuring the word "MILLIPORE" in a blue, sans-serif font above the word "SIGMA" in a larger, bold, blue, sans-serif font.