User Guide

Human High Sensitivity T Cell Panel Serum-Free

96-Well Plate Assay

HSTCSF-28SK HSTCSF28SPMX13 HSTCSF28SPMX21

Introduction2
Principle4
Storage Conditions Upon Receipt5
Reagents Supplied
Materials Required
Safety Precautions
Unique Features of the Human High Sensitivity T Cell Panel Serum-Free 11
Technical Guidelines
Sample Collection and Storage 14 Preparation of Tissue Culture Supernatant14
Preparation of Reagents for Immunoassay14

Preparation of Antibody- Immobilized Beads Preparation of Quality Controls. Preparation of Wash Buffer Preparation of Human High Sensitivity T Cell Standard	14 15
Immunoassay Procedure	18
Plate Washing Solid Plate Filter Plate (Catalogue No. MX-PLATE)	20
Equipment Settings	21
Quality Controls	23
Assay Characteristics Cross-Reactivity	
Troubleshooting	24
Product Ordering	29
Well Map	31
Notice	32
Technical Assistance	32 32



1

Introduction

Cytokines are immunomodulatory polypeptides that play key roles in both adaptive and innate immune responses. A generic term, "cytokines" includes myokines (produced by mononuclear phagocytic cells), lymphokines (produced by activated Th cells), interleukins (acting as mediators between T cells) and chemokines (responsible for T-cell migration). One of the regulatory mechanisms of the immune system, cytokines act at the recognition, activation, or effector phases of an immune response, modulating the development and functional activities of the subtypes of T cells, B cells and myeloid cells. Consequently, research involving cytokines plays a significant role in achieving a deeper understanding of the immune system and its multi-faceted response to most antigens, especially those responses that make up the inflammatory process.

Low levels of inflammation are involved in many clinical and sub-clinical disease states, such as autoimmune disease, cardiovascular disease, diabetes, neurological disorders and cancer. Measuring picogram levels of cytokines, therefore, is critical for understanding the pathogenesis of these diseases.

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
- Detection antibody cocktails designed to yield consistent analyte profiles within panel

In addition, each panel and kit meets stringent manufacturing criteria to ensure batch-to-batch reproducibility. The MILLIPLEX® Human High Sensitivity T Cell Panel Serum-Free thus enables you to focus on the therapeutic potential of cytokines as well as the modulation of even low levels of cytokine expression. 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.

MILLIPLEX® Human High Sensitivity T Cell Panel Serum-Free Magnetic Bead Panel is part of the most versatile system available for cytokines 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 13-plex or 21-plex pre-mixed kit.
- The ability to choose any combination of analytes from our panel of 21 analytes to design a custom 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.

In addition, data obtained from the High Sensitivity T Cell Panel Serum-Free will correlate with data for the respective cytokines in the Human Cytokine/Chemokine Panels I, II and III, furthering your ability to measure specific cytokine response in both normal and disease states.

MILLIPLEX® Human High Sensitivity T Cell Panel Serum-Free is to be used for the simultaneous quantification of any or all of the following analytes in cell/tissue culture supernatant samples: Fractalkine, GM-CSF, IFN γ , IL-1 β , IL-2, IL-4, IL-5, IL-6, IL-7, IL-8, IL-10, IL 12 (p70), IL-13, IL-17A, IL-21, IL-23, ITAC, MIP-1 α , MIP-1 α , MIP-3 α and TNF α .

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

<code>MILLIPLEX®</code> kits are based on the Luminex® <code>xMAP®</code> 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 MILLIPLEX® Analyst 5.1, 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 ≤ -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.
Human High Sensitivity T Cell Standard	1 vial	Lyophilized	HSTC-8028
Human High Sensitivity T Cell Quality Controls 1 and 2	1 vial each	Lyophilized	HSTC-6028
Set of one 96-Well Plate with 2 sealers	1 plate 2 sealers	-	-
Assay Buffer	1 bottle	15 mL	L-ABIR
10X Wash Buffer Note: Contains 0.05% ProClin™	1 bottle	60 mL	L-WB
Human High Sensitivity T Cell Detection Antibodies	1 bottle	5.5 mL	HSTC-1028
Streptavidin-Phycoerythrin	1 bottle	5.5 mL	MC-SAPE7
Bead Diluent	1 bottle	3.5 mL	LBD
Mixing Bottle (not provided with premixed panel)	1 bottle	-	-

Human High Sensitivity T Cell Panel Serum-Free Antibody-Immobilized Premixed Magnetic Beads

Premixed 13-plex Beads	3.5 mL	1 bottle	HSCYPMX13-MAG
Premixed 21-plex Beads	3.5 mL	1 hottle	HSTCPMX21-MAG

Included Human High Sensitivity T Cell Panel Serum-Free Antibody-Immobilized Beads are dependent on customizable selection of analytes within the panel (see next page).

Human High Sensitivity T Cell Panel Serum-Free Antibody-Immobilized Magnetic Beads

			Magnetic Magnetic	Magnetic Magnetic Magneti			(50X concentration, 90 μL	
Bead/Analyte Name	Bead Region	Premixed Beads	Premixed Beads	Available	Catalogue No.			
Anti-Human ITAC Beads	19		~	~	HITAC-MAG			
Anti-Human GM-CSF Beads	20	✓	~	~	HGMCSF-MAG			
Anti-Human Fractalkine Beads	21		~	~	HFKN-MAG			
Anti-Human IFNy Beads	25	~	~	~	HCYIFNG-MAG			
Anti-Human IL-10 Beads	27	~	~	~	HCYIL10-MAG			
Anti-Human MIP-3a Beads	28		~	~	HMIP3A-MAG			
Anti-Human IL-12 (p70) Beads	33	~	~	~	HIL12P70-MAG			
Anti-Human IL-13 Beads	35	~	~	~	HIL13-MAG			
Anti-Human IL-17A Beads	39		~	~	HIL17-MAG			
Anti-Human IL-1β Bead	46	~	~	~	HCYIL1B-MAG			
Anti-Human IL-2 Beads	48	~	~	~	HIL2-MAG			
Anti-Human IL-21 Beads	52		~	~	HIL21-MAG			
Anti-Human IL-4 Beads	53	~	~	~	HIL4-MAG			
Anti-Human IL-23 Beads	54		~	~	HIL23-MAG			
Anti-Human IL-5 Beads	55	~	~	~	HIL5-MAG			
Anti-Human IL-6 Beads	57	~	~	~	HCYIL6-MAG			
Anti-Human IL-7 Beads	61	~	~	~	HIL7-MAG			
Anti-Human IL-8 Beads	63	~	~	~	HCYIL8-MAG			
Anti-Human MIP-1a	72		~	~	HMIP1A-MAG			
Anti-Human MIP-1β	73		~	~	HMIP1B-MAG			
Anti-Human TNFa Beads	75	~	~	~	HCYTNFA-MAG			

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[™], HTS, 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, Cat. Nos. 40-094, 40-095, 40-096, 40-097 or equivalent) or Handheld Magnetic Separation Block (Catalogue No. 40-285 or equivalent).

Note: If a plate washer or handheld magnetic separation block for magnetic beads is not available, one can use a microtiter filter plate (Catalogue No. MX-PLATE) to run the assay using a Vacuum Filtration Unit (Vacuum Manifold Catalogue No. MSVMHTS00 or equivalent with Vacuum Pump Catalogue No. WP6111560 or equivalent).

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	Cat. No.	Label	Label
Human High Sensitivity T Cell Detect Antibodies	HSTC- 1028	(!)	WARNING. 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.
Human High Sensitivity T Cell Quality Control 1 & 2	HSTC- 6028	₹	DANGER. Harmful if swallowed. Causes serious eye damage. Harmful to aquatic life with long lasting effects. Avoid release to the environment. Wear eye 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.
Human High Sensitivity T Cell Standard	HSTC- 8028	₹	DANGER. Harmful if swallowed. Causes serious eye damage. Harmful to aquatic life with long lasting effects. Avoid release to the environment. Wear eye 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.
Assay Buffer	L-ABIR	(!)	WARNING. 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.
Streptavidin- Phycoerythrin	MC-SAPE7		WARNING. 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	<u>(!)</u>	WARNING. May cause an allergic skin reaction. Wear protective gloves. IF ON SKIN: Wash with plenty of soap and water.

Unique Features of the Human High Sensitivity T Cell Panel Serum-Free

Please read this protocol with care as there are several distinctive steps as summarized below:

- When testing tissue culture or other supernatant, the Quality Control and the Standard Vials should be reconstituted and further diluted in the appropriate control culture medium, which will also be used for the background wells.
- 50 µL background, Standard and Quality Controls are added to their appropriate wells on the assay plate.
- 25 μL Sample and 25 μL Assay Buffer are added to the sample wells resulting in a two-fold sample dilution.
- For Quality Control analysis, analyte concentrations DO NOT NEED to be multiplied by the dilution factor.

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 ≤ -20 °C for 1 month and at ≤ -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 μ L Sheath Fluid PLUS in each well and 75 μ 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 background and for reconstitution of standard curve and controls.
- 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 Tissue Culture Supernatant

- Centrifuge the sample to remove debris and assay immediately or aliquot and store samples at ≤ -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:

- \bullet A maximum of 25 μL per well of diluted or neat tissue culture or other media may be used.
- All samples must be stored in polypropylene tubes. DO NOT STORE SAMPLES IN GLASS.

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 70 μ L from each antibody-bead vial to the Mixing Bottle and bring final volume to 3.5 mL with Bead Diluent. Vortex the mixed beads well. Unused portion may be stored at 2-8 °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 70 μ L from each of the 10 bead vials to the Mixing Bottle. Then add 2.8 mL Bead Diluent.

Preparation of Quality Controls

Reconstitute Quality Control 1 (QC1) and Quality Control 2 (QC2) vials with 250 μL appropriate cell culture media. These are the Stock QC Vials. Invert the Stock Vials several times to mix and vortex. Allow the vials to sit for 5-10 minutes. Label two tubes QC1 and QC2 and add 150 μL sample media to each tube. Remove 50 μL from QC1 or QC2 Stock Vials and add to the 150 μL appropriate cell culture media in the respective QC1 and QC2 tubes and vortex. Use these one to four diluted QCs in the assay. Unused portions may be stored at \le -20 °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 $^{\circ}$ C for up to one month.

Preparation of Human High Sensitivity T Cell Standard

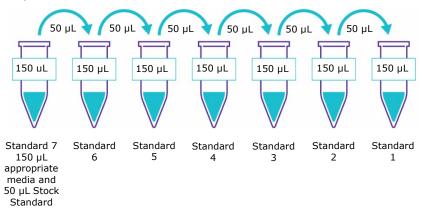
- Prior to use, reconstitute the Human High Sensitivity T Cell Standard with 250 µL appropriate media. Refer to table below for analyte concentrations. Invert the vial several times to mix. Vortex the vial for 10 seconds. Allow the vial to sit for 5-10 minutes. Transfer the reconstituted standard to a polypropylene microfuge tube. This will be used as This is the Stock Standard Vial NOT Standard 7. Unused Standard may be stored at ≤ -20 °C for up to one month.
- 2. Preparation of Working Standards

Label 7 polypropylene microfuge tubes Standard 1 through Standard 7. Add 150 μL of appropriate media to each of the 7 tubes. Prepare serial dilutions by adding 50 μL of the Stock Standard to the Standard 7 tube, mix well and transfer 50 μL of Standard 7 to the Standard 6 tube, mix well and transfer 50 μL of Standard 6 to the Standard 5 tube, mix well and transfer 50 μL of Standard 5 to the Standard 4 tube, mix well and transfer 50 μL of Standard 4 to the Standard 3 tube, mix well and transfer 50 μL of Standard 3 to the Standard 2 tube, mix well and transfer 50 μL of Standard 3 to the Standard 2 tube, mix well and transfer 50 μL of Standard 1 tube and mix well. The 0 pg/mL standard (Background) will be appropriate sample media.

Standard No.	Volume of Appropriate Media to add	Volume of Standard to Add
Stock Standard	250 μL	0

Standard No.	Volume of Appropriate Media to add	Volume of Standard to Add
Standard 7	150 μL	50 μL of Stock Standard
Standard 6	150 μL	50 μL of Standard 7
Standard 5	150 μL	50 μL of Standard 6
Standard 4	150 μL	50 μL of Standard 5
Standard 3	150 μL	50 μL of Standard 4
Standard 2	150 μL	50 μL of Standard 3
Standard 1	150 μL	50 μL of Standard 2

Preparation of Standards



Standard	ITAC, IL-10 (pg/mL)	GM-CSF (pg/mL)	Fractalkine (pg/mL)	IFNy, MIP-3a (pg/mL)	IL-12p70, IL-1β, IL-2, IL-5 (pg/mL)
Standard 1	1.46	1.22	18.3	0.61	0.49
Standard 2	5.86	4.88	73.2	2.44	1.95
Standard 3	23.4	19.5	293.0	9.8	7.81
Standard 4	93.8	78.1	1,171.9	39	31.3
Standard 5	375	312.5	4,687.5	156	125
Standard 6	1,500	1,250	18,750	625	500
Standard 7	6,000	5,000	75,000	2,500	2,000
Standard	IL-13, IL-21 (pg/mL)	IL-17A (pg/mL)	IL-4 (pg/mL)	IL-23 (pg/mL)	IL-6 (pg/mL)
Standard 1	0.24	0.73	1.83	7.93	0.18
Standard 2	0.98	2.93	7.32	31.7	0.73
Standard 3	3.91	11.7	29.3	127.0	2.93
Standard 4	15.63	46.9	117.2	507.8	11.7
Standard 5	62.5	187.5	468.8	2,031.3	46.9
Standard 6	250	750	1,875	8,125	187.5
Standard 7	1,000	3,000	7,500	32,500	750
Standard	IL-7 (pg/mL)	IL-8, MIP-1a (pg/mL)	MIP-1β (pg/mL)	TNFa (pg/mL)	
Standard 1	0.37	0.31	0.92	0.43	
Standard 2	1.46	1.22	3.66	1.71	
Standard 3	5.86	4.88	14.7	6.84	
Standard 4	23.4	19.5	58.6	27.3	
Standard 5	93.8	78.1	234.4	109.4	
Standard 6	375	312.5	937.5	437.5	
Standard 7	1,500	1,250	3,750	1,750	

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 assav.
- 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.

- If using a filter plate, set the filter plate on a plate holder at all times during reagent dispensing and incubation steps so that the bottom of the plate does not touch any surface.
 - Add 200 µL 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).
 - Decant Wash Buffer and remove the residual amount from all wells by inverting the plate and tapping it smartly onto absorbent towels several times.
 - Add 50 µL of each diluted Standard or Quality Control into the appropriate wells (NOT from Stock Vials). The appropriate control culture medium should be used for 0 pg/mL standard (Background).
 - Add 25 μL of Assay Buffer to the sample wells.
 - 5. Add 25 µL of Sample into the sample wells.
 - Vortex Mixing Bottle and add 25 µL of the Mixed or Premixed Beads to each well.

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

 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. Add 200 µL 1X Wash Buffer per well



Shake 10 min, RT Decant

- Add 50 µL Standard or Control to appropriate wells
- Add 50 µL appropriate control culture media to background wells
- Add 25 µL Assay Buffer to sample wells
- Add 25 µL Samples to sample wells
- Add 25 μL Beads to each well



Incubate overnight (16-18 hours) at 2-8 °C

- Gently remove well contents and wash plate 3 times following instructions listed in the **PLATE WASHING** section.
- Add 50 μL of Detection Antibodies into each well.

Note: Allow the Detection Antibodies to warm to room temperature prior to addition.

- Seal, cover with foil and incubate with agitation on a plate shaker for 1 hour at room temperature (20-25 °C). DO NOT ASPIRATE AFTER INCUBATION.
- Add 50 μL Streptavidin-Phycoerythrin to each well containing the 50 μL of Detection Antibodies.
- Seal, cover with foil and incubate with agitation on a plate shaker for 30 minutes at room temperature (20-25 °C).
- Gently remove well contents and wash plate 3 times following instructions listed in the **PLATE WASHING** section.
- 14. Add 150 μ L of Sheath Fluid PLUS (or Drive Fluid PLUS if using MAGPIX®) to all wells. Resuspend the beads on a plate shaker for 5 minutes.
- Run plate on Luminex® 200™, HTS, FLEXMAP 3D®, MAGPIX® with xPONENT® software or xMAP® INTELLIFLEX with INTELLIFLEX Software.
- Save and analyze the Median Fluorescent Intensity (MFI) data using a 5-parameter logistic or spline curvefitting method for calculating analyte concentrations in samples.

Note: Because of the built-in two-fold sample dilution, for all neat samples, multiply the calculated concentrations by two. For two-fold diluted samples, multiply the calculated concentrations by four. Calculated Quality Control concentrations do not require multiplication by a dilution factor.

Remove well contents and wash 3X with 200 µL Wash Buffer

Add 50 µL Detection Antibodies per well



Incubate 1 hour at RT

Do Not Aspirate

Add 50 µL Streptavidin-Phycoerythrin per well



Incubate
30 minutes at RT

Remove well contents and wash 3X with 200 µL Wash Buffer

Add 150 μ L Sheath Fluid PLUS or Drive Fluid PLUS per well

Read on Luminex® instrument (100 µL, 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 μL 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 µL 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.

Filter Plate (Catalogue No. MX-PLATE)

If using a filter plate, use a vacuum filtration manifold to remove well contents. Wash plate with 200 μ L/well of Wash Buffer, removing Wash Buffer by vacuum filtration after each wash. Repeat wash steps as recommended in the Assay Procedure.

Equipment Settings

Luminex® 200™, HTS, 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™ and HTS	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 $^{\otimes}$ 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 μL	
Gate Settings	8,000 to 15,000	
Reporter Gain	Default (low PMT)	
Time Out	60 seconds	
Bead Set	Customizable 21-ple	x Beads
_	ITAC	19
	GM-CSF	20
	Fractalkine	21
	IFNγ	25
	IL-10	27
	MIP-3a	28
	IL-12 (p70)	33
	IL-13	35
	IL-17A	39
	IL-1β	46
	IL-2	48
	IL-21	52
	IL-4	53
	IL-23	54
	IL-5	55
	IL-6	57
	IL-7	61
	IL-8	63
	MIP-1a	72
	MIP-1β	73

TNFa

75

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 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.

Troubleshooting

Problem	Probable Cause	Solution
	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	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.
		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®.

Problem	Probable Cause	Solution					
	Background wells were contaminated	Avoid cross-well contamination by using sealer appropriately and pipetting with multichannel pipettes without touching reagent in plate.					
Background is too high	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.					
	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.					
Beads not	Wrong bead regions in protocol template	Check kit protocol for correct bead regions or analyte selection.					
in region or gate	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.					
Signal for whole plate	Incorrect or no Detection Antibody was added	Add appropriate Detection Antibody and continue.					
is same as background	Streptavidin- Phycoerythrin was not added	Add Streptavidin-Phycoerythrin according to protocol. If Detection Antibody has already been removed, sensitivity may be low.					

Problem Probable Cause		Solution					
Low signal for standard	Detection Antibody may have been removed prior to adding Streptavidin- Phycoerythrin	May need to repeat assay if desired sensitivity not achieved.					
curve	Incubations done at inappropriate temperatures, timings, or agitation	Assay conditions need to be checked.					
Signals too high, standard curves are	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.					
saturated	Plate incubation was too long with standard curve and samples	Use shorter incubation time.					
	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.					
Sample readings are out of range	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.					

Problem	Probable Cause	Solution					
	Multichannel pipette may not be calibrated	Calibrate pipettes.					
	Plate washing was not uniform	Confirm all reagents are removed completely in all wash steps.					
High variation	Samples may have high particulate matter or other interfering substances	See above.					
in samples and/or standards	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.					

For Filter Plates Only

	Vacuum pressure is insufficient	Increase vacuum pressure such that 0.2 mL buffer can be suctioned in 3-5 seconds.					
Filter plate will not vacuum	Samples have insoluble particles	Centrifuge samples just prior to assay set-up and use supernatant.					
	High lipid concentration	After centrifugation, remove lipid layer and use supernatant.					
	Vacuum pressure too high	Adjust vacuum pressure such that 0.2 mL buffer can be suctioned in 3-5 seconds. May need to transfer contents to a new (blocked) plate and continue.					
	Plate set directly on table or absorbent towels during incubations or reagent additions	Set plate on plate holder or raised edge so bottom of filter is not touching any surface.					
Plate leaked	Insufficient blotting of filter plate bottom causing wicking	Blot the bottom of the filter plate well with absorbent towels after each wash step.					
	Pipette touching plate filter during additions	Pipette to the side of plate.					
	Probe height not adjusted correctly	Adjust probe to 3 alignment discs in well H6.					
	Sample too viscous	May need to dilute sample.					

Product Ordering

Order products online at SigmaAldrich.com.

Replacement Reagents	Catalogue Number			
Human High Sensitivity T Cell Standard	HSTC-8028			
Human High Sensitivity T Cell Quality Controls 1 and 2	HSTC-6028			
Human High Sensitivity T Cell Detection Antibodies	HSTC-1028			
Streptavidin-Phycoerythrin	MC-SAPE7			
Assay Buffer	L-ABIR			
Set of two 96-Well plates with sealers	MAG-PLATE			
Bead Diluent	LBD			
10X Wash Buffer	L-WB			

Antibody-Immobilized Magnetic Beads

Bead/Analyte Name	Bead No.	Catalogue Number
ITAC	19	HITAC-MAG
GM-CSF	20	HGMCSF-MAG
Fractalkine	21	HFKN-MAG
IFNγ	25	HCYIFNG-MAG
IL-10	27	HCYIL10-MAG
MIP-3a	28	HMIP3A-MAG
IL-12 (p70)	33	HIL12P70-MAG
IL-13	35	HIL13-MAG
IL-17A	39	HIL17-MAG
IL-1β	46	HCYIL1B-MAG
IL-2	48	HIL2-MAG
IL-21	52	HIL21-MAG
IL-4	53	HIL4-MAG
IL-23	54	HIL23-MAG
IL-5	55	HIL5-MAG
IL-6	57	HCYIL6-MAG
IL-7	61	HIL7-MAG
IL-8	63	HCYIL8-MAG
MIP-1a	72	HMIP1A-MAG
MIP-1β	73	HMIP1B-MAG
TNFa	75	HCYTNFA-MAG
Premixed 13-plex Beads		HSCYPMX13-MAG
Premixed 21-plex Beads		HSTCPMX21-MAG

Well Map

	1	2	3	4	5	6	7	8	9	10	11	12
Α	0 Standard (Background)	Standard 4	QC-1 Control	Etc.								
В	0 Standard (Background	Standard 4	QC-1 Control									
С	Standard 1	Standard 5	QC-2 Control									
D	Standard 1	Standard 5	QC-2 Control									
Е	Standard 2	Standard 6	Sample 1									
F	Standard 2	Standard 6	Sample 1									
G	Standard 3	Standard 7	Sample 2									
Н	Standard 3	Standard 7	Sample 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.

Standard Warranty

The applicable warranty for the products listed in this publication may be found at SigmaAldrich.com/terms.

Contact Information

For the location of the office nearest you, go to SigmaAldrich.com/offices.

Safety Data Sheets (SDS)

Safety Data Sheets are available on the product page at 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.

© 2023 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

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

