

User Guide

# Rat/Mouse GIP (Total) ELISA Kit

## 96-Well Plate

### EZRMGIP-55K

Intended Use.....	2	Calculations .....	12
Principles of Assay.....	2	Interpretation.....	12
Reagents Supplied.....	3	Acceptance Criteria.....	12
Storage and Stability .....	4	Standard Curve .....	13
Reagent Precautions .....	4	Assay Characteristics.....	14
Sodium Azide.....	4	Sensitivity.....	14
Hydrochloric Acid .....	4	Specificity .....	14
Symbol Definitions.....	5	Precision .....	14
Materials Required.....	6	Spike Recovery of Rat GIP in Rat Serum Samples.....	15
Sample Collection and Storage.....	6	Linearity of Sample Dilution .....	16
Sample Preparation .....	7	Quality Controls.....	17
Rat/Mouse GIP		Troubleshooting.....	17
Standard Preparation .....	7	Product Ordering.....	18
Rat/Mouse GIP Quality Controls		Replacement Reagents.....	18
1 and 2 Preparation .....	8	Notice .....	19
Assay Procedure .....	8	Technical Assistance .....	19
Assay Procedure for Rat/Mouse		Terms and Conditions of Sale.....	19
GIP (Total) ELISA Kit .....	10	Contact Information.....	19
Microtiter Plate Arrangement.....	11		

---

## Intended Use

This Rat/Mouse GIP (Total) ELISA kit is used for the non-radioactive quantification of Rat/Mouse GIP (Total) in rat or mouse serum, plasma, tissue extracts or cell culture media samples. This kit has 100% cross reactivity to rat GIP and mouse GIP and measures both GIP(1-42) and GIP(3-42). One kit is sufficient to measure 39 unknown samples in duplicate.

**This kit is for Research Use Only. Not for Use in Diagnostic Procedures.**

## Principles of Assay

This assay is a Sandwich ELISA based, sequentially, on:

- Capture of rat or mouse GIP molecules from samples to the wells of a microtiter plate coated by a pre-titered amount of anti-GIP monoclonal antibodies
- Wash away of unbound materials from samples
- Binding of a second biotinylated anti-GIP polyclonal antibody to the captured molecules
- Wash away of unbound materials from samples
- Incubation of streptavidin-Horseradish peroxidase conjugate to bind to the immobilized biotinylated antibodies
- Wash away of free enzyme conjugates, and
- Quantification of immobilized antibody-enzyme conjugates by monitoring horseradish peroxidase activities in the presence of the substrate 3,3',5,5'-tetramethylbenzidine

The enzyme activity is measured spectrophotometrically by the increased absorbency at 450 nm, corrected from the absorbency at 590 nm, after acidification of formed products. Since the increase in absorbency is directly proportional to the amount of captured rat or mouse GIP in the unknown sample, the latter can be derived by interpolation from a reference curve generated in the same assay with reference standards of known concentrations of Rat GIP.

## Reagents Supplied

Each kit is sufficient to run one 96-well plate and contains the following reagents:

**Note:** Store all reagents at 2-8 °C.

Reagents Supplied	Volume	Quantity	Cat. No.
GIP ELISA Plate with 2 plate sealers <b>Note:</b> Unused strips should be resealed in the foil pouch with the desiccant provided and stored at 2-8 °C.	-	1 plate 2 plates	EP54
10X HRP Wash Buffer Concentrate	50 mL	2 bottles	EWB-HRP
Rat/Mouse GIP Standard	0.5 mL/vial Lyophilized	1 vial	E8055-K
Rat/Mouse GIP Quality Controls 1 and 2	0.5 mL/vial Lyophilized	2 vials	E6055-K
Assay Buffer	12 mL	1 bottle	EABGLP
GIP Detection Antibody	11 mL	1 bottle	E1054
Enzyme Solution	12 mL	1 bottle	EHRP-4
Stop Solution 0.3 M HCl ( <b>Caution:</b> Corrosive solution)	12 mL	1 bottle	ET-TMB
Substrate Solution (Light sensitive, avoid unnecessary exposure to light)	12 mL	1 bottle	ESS-TMB2
Matrix Solutions	1 mL	1 vial	EMTX

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

---

## Storage and Stability

Recommended storage for kit components is 2-8 °C.

All components are shipped and stored at 2-8 °C. Reconstituted standards and controls can be frozen for future use but repeated freeze/thaw cycles should be avoided. Refer to expiration dates on all reagents prior to use. Do not mix reagents from different kits unless they have the same lot numbers.

## Reagent Precautions

### Sodium Azide





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.

### Hydrochloric Acid

Hydrochloric acid is corrosive and can cause eye and skin burns. It is harmful if swallowed and can cause respiratory and digestive tract burns. Avoid contact with skin and eyes. Do not swallow or ingest.

**Note:** See Full Labels of Hazardous Components on next page.

## Symbol Definitions

Ingredient	Cat. No.	Full Label
Rat/Mouse GIP Quality Controls 1 & 2	E6055-K	 <p><b>Warning.</b> Harmful if swallowed. Toxic to aquatic life with long lasting effects. Avoid release to the environment.</p>
Rat/Mouse GIP Standard	E8055-K	 <p><b>Warning.</b> Harmful if swallowed. Toxic to aquatic life with long lasting effects. Avoid release to the environment.</p>
Stop Solution	ET-TMB	 <p><b>Warning.</b> May be corrosive to metals.</p>
10X HRP Wash Buffer Concentrate	EWB-HRP	 <p><b>Warning.</b> May cause an allergic skin reaction. Wear protective gloves. IF ON SKIN: Wash with plenty of soap and water.</p>

---

## Materials Required (Not Provided)

- Multi-channel Pipettes and pipette tips: 5  $\mu$ L-50  $\mu$ L and 50  $\mu$ L-300  $\mu$ L
- Pipettes with tips: 10  $\mu$ L-20  $\mu$ L or 20  $\mu$ L-100  $\mu$ L
- Buffer and Reagent Reservoirs
- Vortex Mixer
- De-ionized water
- Microtiter Plate Reader capable of reading absorbency at 450 nm
- Orbital Microtiter Plate Shaker
- Absorbent Paper or Cloth

## Sample Collection and Storage

**Note:** Although DPPIV inhibitor is not required to be added to serum/plasma samples for measurement of total GIP, we recommend that DPPIV inhibitor be added to the serum/plasma samples during the sample collection so that the same samples could be used in the future for the measurement of intact (1-42) GIP with an assay that is capable of selectively measuring only the intact GIP.

1. To prepare serum samples, whole blood is directly drawn into a centrifuge tube that contains no anti-coagulant. Let blood clot at room temperature for 30 min. Promptly centrifuge the clotted blood at 2,000 to 3,000  $\times g$  for 15 minutes at  $4 \pm 2$  °C. Transfer and store serum samples in separate tubes. Date and identify each sample. Use freshly prepared serum or aliquot and store samples at  $\leq -20$  °C for later use. For long-term storage, keep at  $-70$  °C. Avoid freeze/thaw cycles.
2. To prepare plasma samples, whole blood should be collected into centrifuge tubes containing enough K3EDTA to achieve a final concentration of 1.735 mg/mL and centrifuged immediately after collection. Observe the same precautions in the preparation of serum samples.
3. If heparin is to be used as an anticoagulant, the effect on the assay outcome at the dose of heparin used should be pre-determined.
4. Avoid using samples with gross hemolysis or lipemia.

## Sample Preparation

1. No dilution or preparation is needed for normal serum or plasma samples. In the event that any sample is above 2000 pg/mL range, dilutions should be performed using the Matrix Solution provided.
2. Tissue extracts or cell culture samples may require dilution. Dilutions should be performed using the assay buffer provided.

## Rat/Mouse GIP Standard Preparation

1. Use care in opening the lyophilized Standard vial. Using a pipette, reconstitute the Rat/Mouse GIP Standard with 0.5 mL distilled or deionized water into the vial to give a 2000 pg/mL concentration of Standard. Invert and mix gently, let sit for 5 minutes then mix well.
2. Label five tubes 666.7, 222.2, 74.1, 24.7, and 8.2 pg/mL. Add 100  $\mu$ L Assay Buffer to each of the five tubes. Prepare 3 times serial dilutions by adding 50  $\mu$ L of the 2000 pg/mL reconstituted standard to the 666.7 pg/mL tube, mix well and transfer 50  $\mu$ L of the 666.7 pg/mL standard to the 222.2 pg/mL tube, mix well and transfer 50  $\mu$ L of the 222.2 pg/mL standard to the 74.1 pg/mL tube, mix well and transfer 50  $\mu$ L of the 74.1 pg/mL standard to the 24.7 pg/mL tube, mix well and transfer 50  $\mu$ L of the 24.7 pg/mL standard to the 8.2 pg/mL tube and mix well.

**Note:** Do not use a Repeater pipette. Change tip for every dilution. Wet tip with Standard before dispensing. Unused portions of reconstituted standard should be stored at  $\leq -20$  °C. Avoid multiple freeze/thaw cycles.

<b>Standard Concentration (pg/mL)</b>	<b>Volume of Deionized Water to Add</b>	<b>Volume of Standard to Add</b>
2000	0.5 mL	0

<b>Standard Concentration (pg/mL)</b>	<b>Volume of Assay Buffer to Add</b>	<b>Volume of Standard to Add</b>
666.7	100 $\mu$ L	50 $\mu$ L of 2000 pg/mL
222.2	100 $\mu$ L	50 $\mu$ L of 666.7 pg/mL
74.1	100 $\mu$ L	50 $\mu$ L of 222.2 pg/mL
24.7	100 $\mu$ L	50 $\mu$ L of 74.1 pg/mL
8.2	100 $\mu$ L	50 $\mu$ L of 24.7 pg/mL

## Rat/Mouse GIP Quality Controls 1 and 2 Preparation

Use care in opening the lyophilized Quality Control vials. Using a pipette, reconstitute each of the Rat/Mouse GIP Quality Control 1 and Quality Control 2 with 0.5 mL distilled or deionized water into the vials. Invert and mix gently, let sit for 5 minutes then mix well.

## Assay Procedure

Pre-warm all reagents to room temperature immediately before setting up the assay.

1. Dilute the 10X Wash Buffer 10-fold by adding the entire contents of both bottles of buffer to 900 mL de-ionized or distilled water.
2. Remove required number of strips from the Microtiter Assay Plate. Unused strips should be resealed in the foil pouch with the desiccant provided and stored at 2–8 °C. Assemble strips in an empty plate holder and add 300  $\mu$ L of diluted Wash Buffer to each well. Incubate at room temperature for 5 minutes. Decant wash buffer and remove the residual amount from all wells by inverting the plate and tapping it smartly onto absorbent towels several times. Do not let wells dry before proceeding to the next step. If an automated machine is used for the assay, use a gentle wash program for all washing steps described in this protocol.
3. Add in duplicate, 90  $\mu$ L Assay Buffer to the blank wells and sample wells. (See [Microtiter Plate Arrangement](#) for suggested well orientation).
4. Add in duplicate, 80  $\mu$ L Assay Buffer to Standard wells, QC1 and QC2 wells.
5. Add 10  $\mu$ L Matrix Solution to Blank wells, Standard wells, and QC1 and QC2 wells.
6. Add in duplicate, 10  $\mu$ L Rat/Mouse GIP Standards in the order of ascending concentration to the appropriate wells. Add in duplicate, 10  $\mu$ L QC1 and 10  $\mu$ L QC2 to the appropriate wells. Add sequentially, 10  $\mu$ L of the unknown samples in duplicate to the remaining wells. For best result all additions should be completed within 30 minutes.
7. Cover the plate with plate sealer and incubate at room temperature for 1.5 hours on an orbital microtiter plate shaker set to rotate at moderate speed, approximately 400 to 500 rpm.
8. Remove plate sealer and decant solutions from the plate. Tap as before to remove residual solutions in the wells.
9. Wash wells 3 times with diluted Wash Buffer, 300  $\mu$ L per well per wash. Decant and tap firmly after each wash to remove residual buffer.
10. Add 100  $\mu$ L Detection Antibody to all wells. Cover the plate with plate sealer and incubate at room temperature for 1 hour on an orbital microtiter plate shaker set to rotate at moderate speed, approximately 400 to 500 rpm.
11. Remove plate sealer and decant solutions from the plate. Tap as before to remove residual solutions in the wells.



- 
12. Wash wells 3 times with diluted Wash Buffer, 300  $\mu$ L per well per wash. Decant and tap firmly after each wash to remove residual buffer.
  13. Add 100  $\mu$ L Enzyme Solution to each well. Cover plate with sealer and incubate with moderate shaking at room temperature for 30 minutes on the microtiter plate shaker.
  14. Remove sealer, decant solutions from the plate, and tap plate to remove the residual fluid.
  15. Wash wells 3 times with diluted Wash Buffer, 300  $\mu$ L per well per wash. Decant and tap firmly after each wash to remove residual buffer.
  16. Add 100  $\mu$ L of Substrate Solution to each well, cover plate with sealer and shake on the plate shaker for approximately 5 to 20 minutes (A longer development time may be needed if using a plate washer). Blue color should be formed in wells of GIP standards with intensity proportional to increasing concentrations of GIP.

**Note:** Please be aware that the color may develop more quickly or more slowly than the recommended incubation time depending on the localized room temperature. Please visually monitor the color development to optimize the incubation time. One can monitor color development using 370 nm filter, if available on the spectrophotometer. When the absorbance is between 1.2 and 1.8 at 370 nm, the stop solution can be added to terminate the color development.

17. Remove sealer and add 100  $\mu$ L Stop Solution (**Caution:** Corrosive Solution) and shake plate by hand to ensure complete mixing of solution in all wells. The blue color should turn to yellow after acidification. Wipe the bottom of the microtiter plate to remove any residue prior to reading on plate reader. Read absorbance at 450 nm and 590 nm in a plate reader within 5 minutes and ensure that there are no air bubbles in any well. Record the difference of absorbance units. The absorbance of the highest GIP standard should be approximately 2.0-3.2, or not to exceed the capability of the plate reader used.

## Assay Procedure for Rat/Mouse GIP (Total) ELISA Kit

Well #	Step 1	Step 2	Step 3-4	Step 5	Step 6	Step 7-9	Step 10	Step 11-12	Step 13	Step 14-15	Step 16	Step 17	
A1, B1	Dilute both bottles of 10X Wash Buffer with 900 mL Deionized Water.	Wash plate with 300 $\mu$ L Wash Buffer and incubate at room temperature for 5 minutes. Remove residual buffer by tapping smartly on absorbent towels.	Assay Buffer	Matrix	Standards/ Controls/ Samples	Seal, Agitate, Incubate 1.5 hours at Room Temperature. Wash 3X with 300 $\mu$ L Wash Buffer.	Detection Antibody	Seal, Agitate, Incubate 1 hour at Room Temperature. Remove residual buffer by tapping smartly on absorbent towels. Wash 3X with 300 $\mu$ L Wash Buffer.	Enzyme Solution	Seal, Agitate, Incubate 30 minutes at Room Temperature. Wash 3X with 300 $\mu$ L Wash Buffer.	Substrate	Seal, Agitate, Incubate for 5-20 minutes at Room Temperature.	Stop Solution
C1, D1			90 $\mu$ L	10 $\mu$ L	-		100 $\mu$ L						
E1, F1			80 $\mu$ L	10 $\mu$ L	10 $\mu$ L of 8.2 pg/mL Standard		100 $\mu$ L						
G1, H1			80 $\mu$ L	10 $\mu$ L	10 $\mu$ L of 24.7 pg/mL Standard		100 $\mu$ L						
A2, B2			80 $\mu$ L	10 $\mu$ L	10 $\mu$ L of 74.1 pg/mL Standard		100 $\mu$ L						
C2, D2			80 $\mu$ L	10 $\mu$ L	10 $\mu$ L of 222.2 pg/mL Standard		100 $\mu$ L						
E2, F2			80 $\mu$ L	10 $\mu$ L	10 $\mu$ L of 666.7 pg/mL Standard		100 $\mu$ L						
G2, H2			80 $\mu$ L	10 $\mu$ L	10 $\mu$ L of 2000 pg/mL Standard		100 $\mu$ L						
A3, B3			80 $\mu$ L	10 $\mu$ L	10 $\mu$ L of QC 1		100 $\mu$ L						
C3, D3			80 $\mu$ L	10 $\mu$ L	10 $\mu$ L of QC 2		100 $\mu$ L						
E3, F3			90 $\mu$ L	-	10 $\mu$ L of Sample		100 $\mu$ L						
G3, H3			90 $\mu$ L	-	10 $\mu$ L of Sample		100 $\mu$ L						

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

## Microtiter Plate Arrangement

### Rat/Mouse GIP (Total) ELISA

	1	2	3	4	5	6	7	8	9	10	11	12
A	Blank	222.2 pg/mL	QC 2									
B	Blank	222.2 pg/mL	QC 2									
C	8.2 pg/mL	666.7 pg/mL	Sample									
D	8.2 pg/mL	666.7 pg/mL	Sample									
E	24.7 pg/mL	2000 pg/mL	Sample									
F	24.7 pg/mL	2000 pg/mL	Sample									
G	74.1 pg/mL	QC 1	Sample									
H	74.1 pg/mL	QC 1	Etc.									

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

---

## Calculations

The dose-response curve of this assay fits best to a sigmoidal 5-parameter logistic equation. The results of unknown samples can be calculated with any computer program having a 5-parameter logistic function.

**Note:** When sample volumes assayed differ from 10  $\mu\text{L}$ , an appropriate mathematical adjustment must be made to accommodate for the dilution factor (for example, if 5  $\mu\text{L}$  of sample is used, then calculated data must be multiplied by 2). When sample volume assayed is less than 10  $\mu\text{L}$ , compensate the volume deficit with matrix solution.

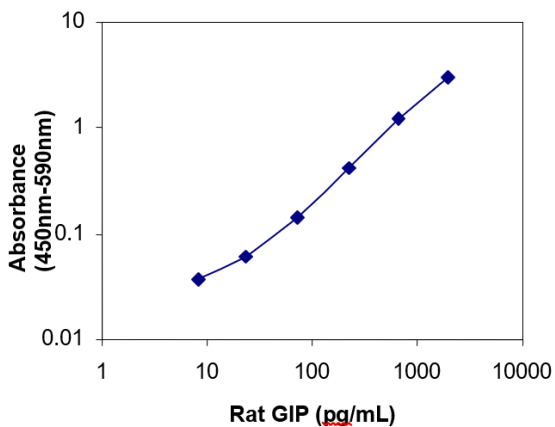
## Interpretation

### Acceptance Criteria

- The assay will be considered accepted when all Quality Control values fall within the calculated Quality Control Range. If any QC's fall outside the control range, review results with a supervisor.
- If the difference between duplicate results of a sample is  $> 10\%$  CV, repeat the sample.
- The limit of sensitivity of this assay is 8.2 pg/mL Rat/Mouse GIP (10  $\mu\text{L}$  sample size).
- The appropriate range of this assay is 8.2 pg/mL to 2000 pg/mL Rat/Mouse GIP (10  $\mu\text{L}$  sample size). Any result greater than 2000 pg/mL in a 10  $\mu\text{L}$  sample should be diluted using matrix solution, and the assay repeated until the results fall within range. Tissue extracts or cell culture media samples greater than 2000 pg/mL in a 10  $\mu\text{L}$  sample should be diluted in Assay Buffer.

## Standard Curve

### Rat/Mouse GIP (Total) ELISA Standard Curve



Typical Standard Curve, not to be used to calculate results.

## Assay Characteristics

### Sensitivity

The lowest level of Rat/Mouse GIP that can be detected by this assay is 8.2 pg/mL when using a 10  $\mu$ L sample size.

### Specificity

The antibody pair used in this assay is specific to rat and mouse GIP and does not significantly cross-react with Glucagon, Oxyntomodulin, GLP-1, and GLP-2.

### Precision

#### Intra-Assay Variation

Sample No.	Mean GIP Levels (pg/mL)	Intra-Assay % CV
1	92	1.0
2	98	2.4
3	173	5.9
4	150	1.1

#### Inter-Assay Variation

Sample No.	Mean GIP Levels (pg/mL)	Inter-Assay % CV
1	80	5.9
2	113	1.1
3	178	4.1

The assay variations of Rat/Mouse GIP ELISA kits were studied on four rat serum samples with varying concentrations of endogenous GIP. The mean intra-assay variation was calculated from results of six duplicate determinations in each assay of the indicated samples. The mean inter-assay variations of each sample were calculated from results of four separate assays with duplicate samples in each assay.

## Spike Recovery of Rat GIP in Rat Serum Samples

Sample	GIP Added (ng/mL)	Expected (ng/mL)	Observed (ng/mL)	% of Recovery
1	0	85	85	100
	49.4	134.4	136	101
	148.9	233.9	238	102
	444.4	529.4	567	107
2	0	93	93	100
	49.4	142.4	146	103
	148.9	241.9	249	103
	444.4	537.4	597	111
3	0	14	14	100
	49.4	63.4	62	98
	148.9	162.9	165	101
	444.4	458.4	512	112

Varying amounts of rat GIP were added to three rat serum samples and the GIP content was determined. The % of recovery = observed GIP concentrations/expected GIP concentrations x 100%.

## Linearity of Sample Dilution

### Effect of Serum Dilution

Sample	Volume Sampled ( $\mu\text{L}$ )	Expected (ng/mL)	Observed (ng/mL)	% of Expected
1	10	43	43	100
	5	22	21	98
	2.5	10.8	11	102
	1.25	5.4	-	-
2	10	62	62	100
	5	31	29	94
	2.5	15.5	14	90
	1.25	7.8	9	116
3	10	21	21	100
	5	10.5	11	105
	2.5	5.25	-	-
	1.25	2.6	-	-

Three rat serum samples with the indicated sample volumes were assayed. Required amounts of matrix were added to compensate for lost volumes below 10  $\mu\text{L}$ . The resulting dilution factors of 1.0, 2.0, 4.0, and 8.0 representing 10  $\mu\text{L}$ , 5  $\mu\text{L}$ , 2.5  $\mu\text{L}$ , and 1.25  $\mu\text{L}$  sample volumes assayed, respectively, were applied in the calculation of observed GIP concentrations. % expected = observed/expected x 100%.



---

## Quality Controls

The ranges for each analyte in Quality Control 1 and 2 are provided on the card insert, or available at our website [SigmaAldrich.com](http://SigmaAldrich.com).

## Troubleshooting

- To obtain reliable and reproducible results the operator should carefully read this manual and fully understand all aspects of each assay step before attempting to run the assay.
- Throughout the assay the operator should adhere strictly to the procedures with good laboratory practice.
- Have all necessary reagents and equipment ready on hand before starting. Once the assay has been started all steps should be completed with precise timing and without interruption.
- Avoid cross contamination of any reagents or samples to be used in the assay.
- Make sure all reagents and samples are added to the bottom of each well.
- Careful and complete mixing of solutions in the well is critical. Poor assay precision will result from incomplete mixing or cross well contamination due to inappropriate mixing.
- Remove any air bubbles formed in the well after acidification of substrate solution because bubbles interfere with spectrophotometric readings.
- Do not let the absorbency reading of the highest standard reach 2.0 units or higher before adding the stop solution.
- High absorbance in the background or blank wells could be due to:
  - Cross well contamination by standard solution or sample.
  - Inadequate washing of the well with Wash Buffer.
  - Overexposure to light after substrate has been added.

## Product Ordering

Products are available for online ordering at [SigmaAldrich.com](https://www.sigmaaldrich.com).

### Replacement Reagents

<b>Reagents</b>	<b>Cat. No.</b>
GIP ELISA Plate	EP54
10X HRP Wash Buffer Concentrate	EWB-HRP
Rat/Mouse GIP Standards	E8055-K
Rat/Mouse GIP Quality Controls 1 and 2	E6055-K
Assay Buffer	EABGLP
Human Leptin Detection Antibody	E1054
Enzyme Solution	EHRP-4
Substrate	ESS-TMB2
Stop Solution	ET-TMB
Matrix Solution	EMTX
10-pack of Rat/Mouse GIP (Total) ELISA kit	EZRMGIP-55BK

---

## 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).

## Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale 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).

---

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.

© 2005-2024 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

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

IFU-EZRMGIP-55K Rev 02/24

20

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.