



www.mesoscale.com®

**Ordering Information**

MSD Customer Service  
 Phone: 1-240-314-2795  
 Fax: 1-301-990-2776  
 Email: CustomerService@mesoscale.com

**Scientific Support**

Phone: 1-240-314-2798  
 Email: ScientificSupport@mesoscale.com

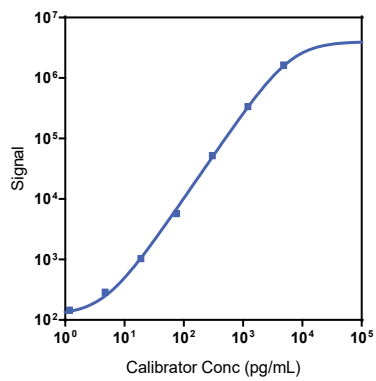
**Company Address**

MESO SCALE DISCOVERY®  
 A division of  
 Meso Scale Diagnostics, LLC.  
 1601 Research Boulevard  
 Rockville, MD 20850-3173 USA

| Product Options     | Catalog Number  | Description   |
|---------------------|---|---|
| <b>Multiplex</b>    | K15067M, K25067M  | U-PLEX Biomarker Group 1 (human)                      |
|                     | K151AEM, K251AEM  | U-PLEX Immuno-Oncology Group 1 (human)                |
|                     | K151ACM, K251ACM  | U-PLEX Metabolic Group 1 (human)                      |
| <b>Singleplex</b>   | K151UDK-1/-2/-4   | U-PLEX Human Eotaxin Assay with SECTOR™ plates        |
|                     | K151UDK-21  | U-PLEX Human Eotaxin Assay with QuickPlex® APT plates |
|                     | K251UDK-2/-4  | U-PLEX Human Eotaxin Assay with 384-well plates       |
| <b>Antibody Set</b> | B21UD-2/-3  | U-PLEX Human Eotaxin Antibody Set                     |
| <b>Protocol</b>     | U-PLEX Product Inserts are available at <a href="http://www.mesoscale.com">http://www.mesoscale.com</a> |   |

The U-PLEX® platform was designed to provide ultimate flexibility for detection of biomarkers in a wide variety of sample types. This datasheet provides the representative performance of the U-PLEX Human Eotaxin Assay tested on U-PLEX 96-well SECTOR plates run as a multiplex. The data do not represent the product specifications. Under your experimental conditions, the assay may perform differently from the representative data. U-PLEX assays are offered in either singleplex or multiplex; both are available on 96- or 384-well plates. See a U-PLEX product insert for instrument compatibility.

**Representative Calibration Curve and Sensitivity**



| Assay   | Median LLOD (pg/mL) | LLOD Range (pg/mL) |
|---------|---------------------|--------------------|
| Eotaxin | 8.2                 | 3.6–11.3           |

The Calibrator curve was fitted with a 4-parameter logistic model with a 1/Y<sup>2</sup> weighting. The lower limit of detection (LLOD) is a calculated concentration corresponding to 2.5 standard deviations above the background (zero Calibrator).

**Precision**

| Control | Average Conc. (pg/mL) | Average Intra-run Conc. (%CV) | Inter-run Conc. (%CV) |
|---------|-----------------------|-------------------------------|-----------------------|
| High    | 961                   | 2.6                           | 7.7                   |
| Mid     | 97                    | 3.1                           | 9.7                   |
| Low     | 9.0                   | 14.7                          | 26.3                  |

Controls were made by spiking Calibrator into assay diluent at 3 levels within the quantitative range of the assay. Average intra-run concentration %CV is the average %CV of the control replicates within an individual run. Inter-run concentration %CV is the variability of controls across multiple runs.

For Research Use Only.  
 Not for use in diagnostic procedures.

# MSD® U-PLEX Human Eotaxin

## Tested Samples

| Sample Type    | Serum (N=10) | Plasma (N=10) | Spiked Plasma (N=5) | Spiked Serum (N=5) |
|----------------|--------------|---------------|---------------------|--------------------|
| Median (pg/mL) | 46           | 101           | 91                  | 52                 |
| Range (pg/mL)  | 12-143       | 26-249        | 40-169              | 49-282             |
| % Detected     | 100          | 100           | 100                 | 100                |

Normal serum and plasma samples were diluted 2-fold prior to testing in the assay.

## Dilution Linearity

| Serum         |                    |                  | EDTA Plasma   |                    |                  |
|---------------|--------------------|------------------|---------------|--------------------|------------------|
| Fold Dilution | Average % Recovery | % Recovery Range | Fold Dilution | Average % Recovery | % Recovery Range |
| 2             | 112                | 107-115          | 2             | 125                | 118-134          |
| 4             | 112                | 108-116          | 4             | 137                | 125-153          |
| 8             | 106                | 100-114          | 8             | 127                | 100-158          |

Normal human serum and EDTA plasma were spiked with Calibrator and tested at different dilutions. Two-fold diluted samples were tested to determine the expected concentration of the analyte. Samples may benefit from additional dilution with assay diluent to reduce matrix effects.

$$\% \text{ Recovery} = (\text{measured concentration} / \text{expected concentration}) \times 100$$

## Spike Recovery

| Spike Level | Serum              |                  | EDTA Plasma        |                  |
|-------------|--------------------|------------------|--------------------|------------------|
|             | Average % Recovery | % Recovery Range | Average % Recovery | % Recovery Range |
| High        | 98                 | 92-101           | 88                 | 80-98            |
| Mid         | 106                | 96-114           | 96                 | 87-105           |
| Low         | 102                | 96-108           | 101                | 95-106           |

Normal serum and plasma were spiked with Calibrator at 3 levels. Two-fold diluted samples were tested to determine the expected concentration of the analyte. Samples may benefit from additional dilution with assay diluent to reduce matrix effects.

$$\% \text{ Recovery} = (\text{measured concentration} / \text{expected concentration}) \times 100$$

## Specificity

The Eotaxin Antibody Set was tested against all of the analytes in Biomarker Group 1, Metabolic Group 1, and Immuno-Oncology Group 1. Any non-specific binding greater than 2.0% is noted below. The U-PLEX Assay Designer shows compatible assays.

$$\% \text{ Nonspecificity} = (\text{nonspecific signal} / \text{specific signal}) \times 100$$

## Diluent Compatibility

Diluents 57 and 3 are provided when this product is ordered in singleplex and when multiplexed with other Biomarker Group 1 assays. Other diluents may be provided when combined with assays from other U-PLEX Groups. See the appropriate Product Insert for details.

## Assay Components

**Calibrator:** Eotaxin is included in Calibrator 2. The Eotaxin Calibrator is a full-length recombinant protein expressed in *E. coli*.

**Antibodies:** The U-PLEX Human Eotaxin Assay uses a mouse monoclonal antibody for capture and a mouse monoclonal antibody for detection.

**Assay generation:** B

**Note:** This datasheet contains representative assay performance data. In custom multiplex formats, the assay may perform differently from the representative data shown.

