



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

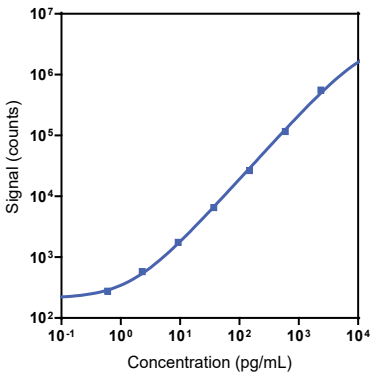
Company Address

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

| Product Options | Catalog Number | Description |
|-----------------|--|---|
| Multiplex | K151AEM, K251AEM | U-PLEX Immuno-Oncology Group 1 (human) |
| Singleplex | K151AQTk-1/-2/-4 | U-PLEX Human proMMP-10 Assay with SECTOR™ plates |
| | K151AQTk-21 | U-PLEX Human proMMP-10 Assay with QuickPlex® plates |
| | K251AQTk-2/-4 | U-PLEX Human proMMP-10 with 384-well plates |
| Antibody Set | B21AQT-2/-3 | U-PLEX Human proMMP-10 Antibody Set |
| Protocol | U-PLEX Product Inserts are available at www.mesoscale.com | |

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 proMMP-10 Assay tested on U-PLEX 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) |
|-----------|---------------------|--------------------|
| proMMP-10 | 0.38 | 0.31–0.55 |

The Calibrator curve was fitted with a 4-parameter logistic model with a 1/Y² 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 | 1,357 | 2.5 | 16.4 |
| Mid | 288 | 3.1 | 12.6 |
| Low | 69 | 1.8 | 13.8 |

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 proMMP-10

Tested Samples

| Sample Type | Serum (N = 8) | EDTA Plasma (N = 8) | Citrate Plasma (N = 8) | Normal Lysate (N = 5) | Tumor Lysate (N = 15) |
|----------------|------------------|------------------------|---------------------------|--------------------------|--------------------------|
| Median (pg/mL) | 272 | 6.0 | 375 | 2.5 | 2.6 |
| Range (pg/mL) | 74–588 | ND–6.0 | 105–819 | 1.1–2.6 | 1.2–3.7 |
| % Detected | 100 | 13 | 100 | 100 | 100 |

Normal serum and plasma samples were diluted 4-fold prior to the assay. Lysates were tested at a protein concentration of 0.5 mg/mL. ND = non-detectable (<LLOD). EDTA and citrate in samples may adversely affect this assay and are not recommended.

Dilution Linearity

| Serum | | | EDTA Plasma | | |
|---------------|--------------------|------------------|---------------|--------------------|------------------|
| Fold Dilution | Average % Recovery | % Recovery Range | Fold Dilution | Average % Recovery | % Recovery Range |
| 2 | 135 | 121–158 | 2 | ND | ND |
| 8 | 90 | 72–108 | 8 | ND | ND |
| 16 | 76 | 67–107 | 16 | ND | ND |

Normal human serum and EDTA plasma were spiked with Calibrator and tested at different dilutions. Percent recovery at each dilution level was normalized to the dilution-adjusted, 4-fold concentration. Samples may benefit from additional dilution with assay diluent to reduce matrix effects. ND = non-detectable (<LLOD).

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

Spike Recovery

| Spike Level | Serum | | EDTA Plasma | |
|-------------|--------------------|------------------|--------------------|------------------|
| | Average % Recovery | % Recovery Range | Average % Recovery | % Recovery Range |
| High | 104 | 82–123 | ND | ND |
| Mid | 102 | 89–116 | ND | ND |
| Low | 104 | 94–115 | ND | ND |

Normal serum and plasma were spiked with Calibrator at 3 levels. Spiked samples were diluted 4-fold to determine the expected concentration of the analyte. Samples may benefit from additional dilution with assay diluent to reduce matrix effects. ND = non-detectable (<LLOD).

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

Specificity

The proMMP-10 Antibody Set was tested for nonspecific binding against all of the analytes in the Immuno-Oncology Group 1 and the majority of analytes in Biomarker Group 1. Any cross-reactivity greater than 2.0% is noted below. The U-PLEX Assay Designer shows compatible assays.

The proMMP-10 assay cross-reacts with the MMP-10 assay as expected. We do not recommend multiplexing these assays on the same plate.

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

Diluent Compatibility

Diluents 58 and 3 are provided when this is ordered in singleplex and multiplex assays.

Assay Components

Calibrator: proMMP-10 is included in Calibrator 32. The human proMMP-10 Calibrator is MMP-10 (18–476) recombinant protein expressed in a mouse cell line.

Antibodies: The U-PLEX Human proMMP-10 Assay uses a mouse monoclonal antibody for capture and a goat polyclonal antibody for detection.

Assay generation: A

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

