



www.mesoscale.com[®]

Ordering Information

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

Scientific Support

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

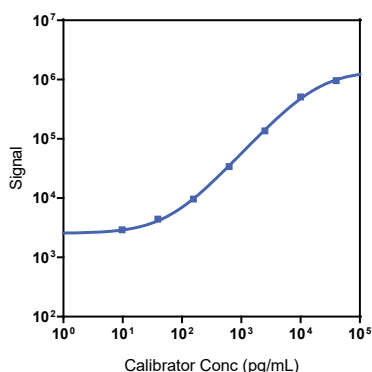
Company Address

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

| Product Options | Catalog Number | Description |
|-----------------|--|--|
| Multiplex | K151ADM, K251ADM | U-PLEX Biomarker Group 2 (human) |
| Singleplex | K151XWK-1/-2/-4 | U-PLEX Human TGF- β 1 Assay with SECTOR [™] plates |
| | K151XWK-21/-22/-24 | U-PLEX Human TGF- β 1 Assay with QuickPlex [®] plates |
| | K251XWK-2/-4 | U-PLEX Human TGF- β 1 Assay with 384-well plates |
| Antibody Set | B20XW-2/-3 | U-PLEX TGF- β 1 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 TGF- β 1 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) |
|----------------|---------------------|--------------------|
| TGF- β 1 | 9.1 | 5.0-10 |

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 | 1,570 | 4.7 | 11.1 |
| Mid | 388 | 5.5 | 11.6 |
| Low | 104 | 5.7 | 16.8 |

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 Assays

Tested Samples

| Sample Type | Serum (N=10) | Plasma (N=10) |
|----------------|--------------|---------------|
| Median (pg/mL) | 1,160 | 2,130 |
| Range (pg/mL) | 426-10,700 | 864-5,110 |
| % Detected | 100 | 100 |

Normal human serum and EDTA plasma samples were tested without dilution prior to the assay. Samples were prepared using an acidification step.

Dilution Linearity

| Serum | | | EDTA Plasma | | | Cell Culture Media | | |
|---------------|--------------------|------------------|---------------|--------------------|------------------|--------------------|--------------------|------------------|
| Fold Dilution | Average % Recovery | % Recovery Range | Fold Dilution | Average % Recovery | % Recovery Range | Fold Dilution | Average % Recovery | % Recovery Range |
| 2 | 136 | 125-157 | 123 | 105-143 | 114 | 105-125 | 136 | 125-157 |
| 4 | 148 | 131-169 | 134 | 108-152 | 114 | 104-125 | 148 | 131-169 |
| 8 | 155 | 142-171 | 139 | 106-154 | 118 | 106-139 | 155 | 142-171 |

Normal human serum, EDTA plasma, and cell culture media were spiked with Calibrator and tested at different dilutions. Undiluted samples were tested to determine the expected concentration of the analyte. Samples may benefit from additional dilution with assay diluent to reduce matrix effects.

% Recovery = (measured concentration / expected concentration) x 100

Spike Recovery

| Spike Level | Serum | | EDTA Plasma | | Cell Culture Media | |
|-------------|--------------------|------------------|--------------------|------------------|--------------------|------------------|
| | Average % Recovery | % Recovery Range | Average % Recovery | % Recovery Range | Average % Recovery | % Recovery Range |
| High | 72 | 61-86 | 57 | 51-67 | 70 | 61-86 |
| Mid | 70 | 64-81 | 58 | 48-68 | 69 | 63-81 |
| Low | 67 | 59-78 | 53 | 47-65 | 66 | 59-78 |

Normal human serum, EDTA plasma, and cell culture media were spiked with Calibrator at 3 levels. Undiluted samples were tested to determine the expected concentration of the analyte. Samples may benefit from additional dilution with assay diluent to reduce matrix effects.

% Recovery = (measured concentration / expected concentration) x 100

Specificity

To assess specificity, the TGF- β 1 Antibody Set was tested individually against a larger panel of analytes for nonspecific binding (TGF- β 1, TGF- β 2, TGF- β 3). Nonspecific binding was less than 0.5%.

% Nonspecificity = (nonspecific signal / specific signal) x 100

It is recommended that acid-treated samples are used for evaluation of TGF- β 1. Samples may benefit from an additional dilution prior to measurement to ensure TGF- β 1 levels are in the quantitative range of the assay.

Diluent Compatibility

Diluents 57 and 3 are provided with this assay. MSD offers a range of assay and antibody diluents for separate purchase. Depending on assay needs, customers may wish to test other diluents.

Assay Components

Calibrator: Human TGF- β 1 is included in Calibrator 11. The TGF- β 1 Calibrator is a full-length recombinant protein expressed in *E. coli*.

Antibodies: The U-PLEX Human TGF- β 1 Assay uses a mouse monoclonal antibody for capture and a chicken polyclonal antibody for detection.

Assay generation: A

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

