



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

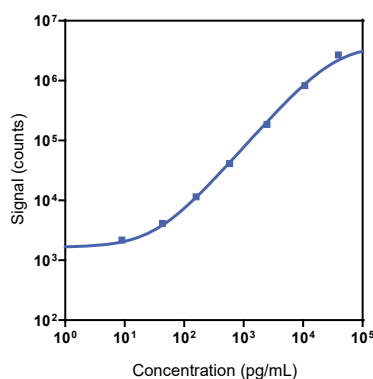
Company Address

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

Product Options	Catalog Number	Description
Multiplex	K152ADM, K252ADM	U-PLEX Biomarker Group 2 (mouse)
Singleplex	K152XUK-1/-2/-4	U-PLEX Mouse TGF- β 2 Assay with SECTOR [™] plates
	K152XUK-21	U-PLEX Mouse TGF- β 2 Assay with QuickPlex [®] APT plates
	K252XUK-2/-4	U-PLEX Mouse TGF- β 2 Assay with 384-well plates
Antibody Set	B20XU-2/-3	U-PLEX Mouse TGF- β 2 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 Mouse TGF- β 2 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 in 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 2	2.5	2.0-2.6

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

Precision

Control	Average Conc. (pg/mL)	Average Intra-run Conc. (%CV)	Inter-run Conc. (%CV)
High	5,140	8.4	11.9
Mid	519	8.6	11.9
Low	72	8.3	11.5

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

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.

MSD® U-PLEX Assays

Tested Samples

Sample Type	Serum (N=13)	Plasma (N=10)
Median (pg/mL)	95	101
Range (pg/mL)	28-218	36-120
% Detected	100	100

Normal serum, EDTA plasma, and cell culture media were diluted 2-fold before 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	130	123-140	2	119	116-122	2	93	87-100
4	139	129-154	4	134	131-137	4	88	83-92
8	147	128-167	8	139	125-146	8	81	75-87

Normal mouse serum and EDTA plasma 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.

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

Spike Recovery

Spike Level	Serum		EDTA Plasma		Cell Culture Media	
	Average % Recovery	% Recovery Range	Average % Recovery	% Recovery Range	Average % Recovery	% Recovery Range
High	53	46-58	56	53-61	117	112-125
Mid	51	47-56	54	51-59	113	109-116
Low	50	47-52	51	49-56	114	109-118

Normal serum and plasma 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.

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

Specificity

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

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

It is recommended that acid-treated samples are used for the evaluation of TGF- β 2. TGF- β 2 assays may experience a modest increase in signal when multiplexed with TGF- β 3 assay due to interaction between capture antibodies. Samples may require additional dilution with assay diluent to reduce matrix effects.

Diluent Compatibility

The data included in this document have been collected with Assay Diluent 41 and Antibody Diluent 45. 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: Mouse TGF- β 2 is included in Calibrator 11. The TGF- β 2 Calibrator is a full-length recombinant protein expressed in *E. coli*.

Antibodies: The U-PLEX Mouse TGF- β 2 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 than the representative data shown.

