



www.mesoscale.com®

Ordering Information

MSD Customer Service
Phone: 1-240-314-2798
: 1-301-990-2776
Email: CustomerService@mesoscale.com

Scientific Support

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

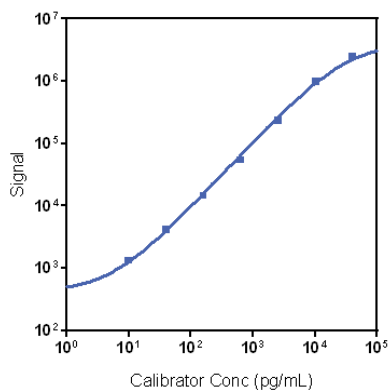
Company Address

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	K151XVK-1/-2/-4	U-PLEX Human TGF-β3 Assay with SECTOR™ plates
	K151XVK-21/-22/-24	U-PLEX Human TGF-β3 Assay with QuickPlex® plates
	K251XVK-2/-4	U-PLEX Human TGF-β3 Assay with 384-well plates
Antibody Set	B20XV-2/-3	U-PLEX TGF-β3 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-β3 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-β3	1.4	1.1-1.4

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	4,010	4.1	102
Mid	425	5.0	107
Low	54	4.8	13.4

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)	ND	ND
Range (pg/mL)	ND	ND
% Detected	0	0

Normal human serum and EDTA plasma samples were diluted 2-fold prior to the assay. Samples were prepared using an acidification step. ND = non-detectable (< LLOD)

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	127	124-132	107	97-121	83	78-88	127	124-132
4	145	137-154	113	98-145	77	73-83	145	137-154
8	154	149-163	113	94-147	71	62-81	154	149-163

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	115	103-119	73	61-89	115	103-119
Mid	108	100-116	74	61-83	108	100-116
Low	105	100-112	72	61-88	105	100-112

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-β3 Antibody Set was tested individually against a larger panel of analytes for nonspecific binding (TGF-β1, TGF-β2, TGF-β3). 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-β3. Samples may benefit from an additional dilution prior to measurement to ensure TGF-β3 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-β3 is included in Calibrator 11. The TGF-β3 Calibrator is a full-length recombinant protein expressed in *E. coli*.

Antibodies: The U-PLEX Human TGF-β3 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.

