

U-PLEX[®] Mouse TGF- β 3 Assay



www.mesoscale.com[®]

Ordering Information

MSD[®] Customer Service
 Phone: 1-301-947-2085
 Fax: 1-301-990-2776
 Email: CustomerService@mesoscale.com

Scientific Support

Phone: 1-301-947-2025
 Email: ScientificSupport@mesoscale.com

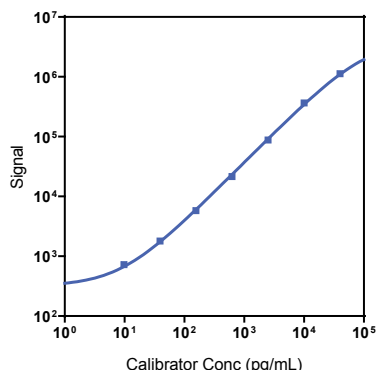
Company Address

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

Product Options	Available in: TGF- β Combo (ms) K15242K
	Individual assay: K152XVK
	Antibody Set: B20XV
Assay Protocol	Refer to the U-PLEX Biomarker Group 2 product insert available at www.mesoscale.com/U-PLEX-documents .

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 3 Assay tested on U-PLEX plates run as a multiplex. The data were generated during the development of the assay and do not represent the product specifications. Under your experimental conditions and with your specific multiplex, the assay may perform differently than the representative data shown. U-PLEX assays are available in multiplex format with other compatible assays. The same assay can also be used to detect a single analyte on MSD GOLD[™] Small Spot Streptavidin plates.

Representative Calibration Curve and Sensitivity



Assay	Median LLOD (pg/mL)	LLOD Range (pg/mL)
TGF- β 3	2.5	1.6-2.9

The calibration curves used to calculate analyte concentrations were established by fitting the signals from the Calibrators using a 4-parameter logistic (or sigmoidal dose-response) model with a $1/Y^2$ weighting. Analyte concentrations were determined from the electrochemiluminescence signals by back-fitting to the calibration curve. The lower limit of detection (LLOD) is a calculated concentration corresponding to the signal 2.5 standard deviations above the background (zero Calibrator).

Precision

	Control	Average Conc. (pg/mL)	Average Intra-run Conc. %CV	Inter-run Conc. %CV
TGF- β 3	High	3,649	8.7	12.6
	Mid	391	9.1	13.1
	Low	53	10.2	14.4

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	Plasma
Median (pg/mL)	16	9
Range (pg/mL)	0-21	0-14
% Detected	67	67

Samples were prepared using an acidification step. No TGF-β3 was detected in samples not treated with acid.

Normal mouse serum and EDTA plasma samples were tested without dilution prior to the assay.

Dilution Linearity

	Serum			Plasma			Cell Culture Media		
	Fold Dilution	Average % Recovery	% Recovery Range	Fold Dilution	Average % Recovery	% Recovery Range	Fold Dilution	Average % Recovery	% Recovery Range
TGF-β3	2	140	137-144	2	130	126-133	2	87	82-92
	4	162	156-168	4	152	143-158	4	80	75-84
	8	177	161-190	8	160	144-169	8	72	65-76

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

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

Spike Recovery

	Spike Level	Serum		Plasma		Cell Culture Media	
		Average % Recovery	% Recovery Range	Average % Recovery	% Recovery Range	Average % Recovery	% Recovery Range
TGF-β3	High	46	40-50	57	53-61	133	131-137
	Mid	46	45-49	57	54-61	130	126-135
	Low	47	44-49	57	54-60	130	127-133

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

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

Specificity

To assess specificity, TGF-β3 Antibody Set was tested individually against a larger panel of recombinant mouse analytes for nonspecific binding (EPO, GM-CSF, IFN-γ, IL-1β, IL-2, IL-4, IL-5, IL-6, IL-9, IL-10, IL-12/IL-23p40, IL-12p70, IL-13, IL-15, IL-16, IL-17A, IL-17A/F, IL-17C, IL-17E/IL-25, IL-17F, IL-21, IL-22, IL-23, IL-27p28/IL-30, IL-31, IL-33, IP-10, KC/GRO, MCP-1, MIP-1α, MIP-1β, MIP-2, MIP-3α, TNF-α, and VEGF-A). Nonspecific binding was less than 0.5%.

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

Diluent Compatibility

The data included in this document has been collected using 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: Human TGF-β3 is included in Calibrator 11 blend. The human TGF-β3 Calibrator is a full-length recombinant human protein expressed in *E. coli*.

Antibodies: The U-PLEX Mouse TGF-β3 Assay uses mouse monoclonal antibody for capture and 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.

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, www.mesoscale.com, MSD, MSD GOLD, MULTI-ARRAY, MULTI-SPOT, SULFO-TAG, U-PLEX, 96 WELL 1, 4, 7, 9, & 10-SPOT (designs), MSD (design), U-PLEX (design), It's All About U, and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.

©2016, 2019 Meso Scale Diagnostics, LLC. All rights reserved.