

U-PLEX[®] Mouse IL-17A/F 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

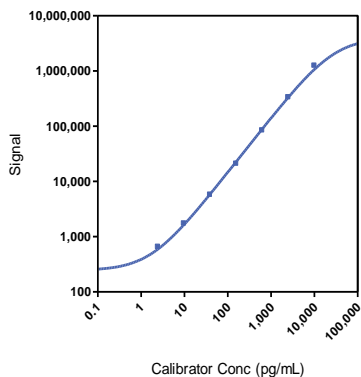
Company Address

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

Product Options	Available as part of U-PLEX Biomarker Group 1 (ms) multiplex combination: K15069L-1/-2/-4
	Individual assay: K152VYK-1/-2/-4; Antibody Set: B22VY-2/B22VY-3
	For more ordering options, please visit www.mesoscale.com
Instrument Compatibility	SECTOR [®] Imager 2400, SECTOR Imager 6000, MESO [®] SECTOR S 600, MESO QuickPlex [®] SQ 120
Sample Type	Mouse serum, EDTA plasma, and cell culture supernatants
Assay Protocol	Refer to the U-PLEX Biomarker Group 1 (Mouse) 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 Mouse IL-17A/F 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)
IL-17A/F	0.61	0.55-0.87

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
IL-17A/F	High	1,590	2.8	7.0
	Mid	212	3.1	11.1
	Low	29	4.3	15.7

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

Spike Recovery

	Spike Level	Serum		Plasma		Cell Culture Media	
		Average % Recovery	% Recovery Range	Average % Recovery	% Recovery Range	Average % Recovery	% Recovery Range
IL-17A/F	High	76	73-80	68	53-74	91	78-107
	Mid	78	69-85	72	57-78	94	76-110
	Low	79	57-95	70	53-77	93	75-109

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

Tested Samples

Sample Type	Serum	Plasma	Stimulated Sample
Median (pg/mL)	ND	3.1	24
Range (pg/mL)	ND-3.2	2.3-4.4	4.0-76
% Detected	25	100	100

ND = non-detectable above the LLOD

Normal mouse serum and EDTA plasma samples were tested without dilution prior to the assay. Spiked serum and spiked plasma represent samples that were spiked with Calibrator and/or cell culture supernatants derived from mouse splenocytes stimulated with different compounds in vitro.

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
IL-17A/F	2	115	109-123	2	123	114-135	2	106	100-114
	4	125	121-129	4	132	123-156	4	98	89-105
	8	125	114-130	8	140	129-166	8	101	91-110

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

Specificity

To assess specificity, IL-17A/F Antibody Set was tested individually against a larger panel of recombinant mouse analytes for nonspecific binding (EPO, GM-CSF, IL-1 β , 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/F, IL-17C, IL-17E, IL-21, IL-22, IL-23, IL-31, IL-33, IP-10, KC/GRO, MCP-1, MIP-1 α , MIP-1 β , MIP-2, MIP-3 α , TGF- β 1, TGF- β 2, TGF- β 3, TNF- α , and VEGF-A). Nonspecific binding was less than 0.5%. IL-17A/F will bind to both IL-17A and IL-17F spots. Do not multiplex with IL-17A. Do not blend Calibrator 7 (IL-17A) with Calibrator 12 (IL-17A/F).

Nonspecificity = (nonspecific signal / specific signal) x 100

Diluent Compatibility

The data included in this document has been collected using Diluents 41 and 45. MSD offers a range of assay and antibody diluents for separate purchase. Depending on your assay needs, other diluents may be tested.

Assay Components

Calibrator: Mouse IL-17A/F is included in Calibrator 12 blend. The full-length recombinant mouse protein expressed in E.coli is used.

Antibodies: The U-PLEX Mouse IL-17A/F Assay uses rat monoclonal antibody for capture and rat monoclonal 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, MSD, MSD GOLD, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, U-PLEX, S-PLEX, V-PLEX, STREPTAVIDIN GOLD, MESO, www.mesoscale.com, SMALL SPOT (design), 96 WELL 1, 4, 7, 9, & 10-SPOT (designs), 384 WELL 1 & 4-SPOT (designs), MSD (design), U-PLEX (design), S-PLEX (design), V-PLEX (design), It's All About U, and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.

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

