

Mouse IL-27p28/IL-30



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

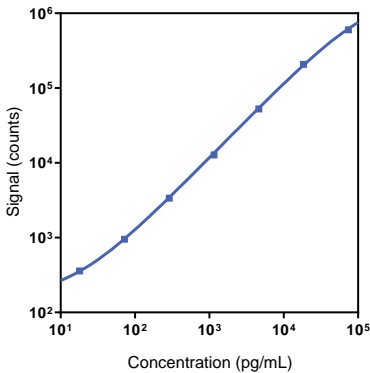
Company Address

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

Product Options	Catalog Number	Description
Multiplex	K15069M, K25069M	U-PLEX Biomarker Group 1 (ms)
	K152ACM, K252ACM	U-PLEX Metabolic Group 1 (ms)
Singleplex	K152WCK-1/-2/-4	U-PLEX Mouse IL-27p28/IL-30 Assay with SECTOR™ plates
	K152WCK-21/-22/-24	U-PLEX Mouse IL-27p28/IL-30 Assay with QuickPlex® plates
	K252WCK-2/-4	U-PLEX Mouse IL-27p28/IL-30 Assay with 384-well plates
Antibody Set	B22WC-2/-3	U-PLEX Mouse IL-27p28/IL-30 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 IL-27p28/IL-30 Assay tested on U-PLEX 96-well 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)
IL-27p28/IL-30	8.7	6.0-11

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 2.5X the standard deviations above the background (zero Calibrator).

Precision

Control	Average Conc. (pg/mL)	Average Intra-run Conc. (%CV)	Inter-run Conc. (%CV)
High	34,100	3.3	3.3
Mid	3,550	2.7	3.2
Low	300	2.2	5.9

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 Mouse IL-27p28/IL-30

Tested Samples

Sample Type	Serum (N=8)	Plasma (N=8)	Stimulated Sample (N=8)
Median (pg/mL)	18	11	22
Range (pg/mL)	13-24	ND-14	ND-993
% Detected	100	63	63

Normal serum and plasma samples were tested without dilution prior to the assay. ND = non-detectable (<LLOD)

Dilution Linearity

Serum			EDTA Plasma		
Fold Dilution	Average % Recovery	% Recovery Range	Fold Dilution	Average % Recovery	% Recovery Range
2	114	108-124	2	115	113-119
4	123	118-131	4	128	124-132
8	127	117-131	8	133	124-140

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.

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

Spike Recovery

Spike Level	Serum		EDTA Plasma	
	Average % Recovery	% Recovery Range	Average % Recovery	% Recovery Range
High	67	65-70	63	61-68
Mid	65	62-66	61	58-63
Low	62	60-64	59	56-63

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.

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

Specificity

To assess specificity, the IL-27p28/IL-30 Antibody Set was tested individually against a larger panel of analytes for nonspecific binding (6CKine/Ccl21, BAFF, BCA-1/BLC, CD40, Eotaxin, EPO, GM-CSF, IFN- α , IFN- β , IFN- γ , 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, 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, MCP-5/Ccl12, MDC, MIP-1 α , MIP-1 β , MIP-2, MIP-3 α , MMP-9 (total), NGAL/LCN2, RANTES, SDF-1 α , TARC, TNF-RI, TNF- α , VEGF-A). Nonspecific binding was less than 0.5%.

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

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 your assay needs, other diluents may be tested.

Assay Components

Calibrator: IL-27p28/IL-30 is included in Calibrator 8. The IL-27p28/IL-30 Calibrator is a full-length recombinant protein expressed in a mouse cell line.

Antibodies: The U-PLEX Mouse IL-27p28/IL-30 Assay uses a goat polyclonal 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 from the representative data shown.

