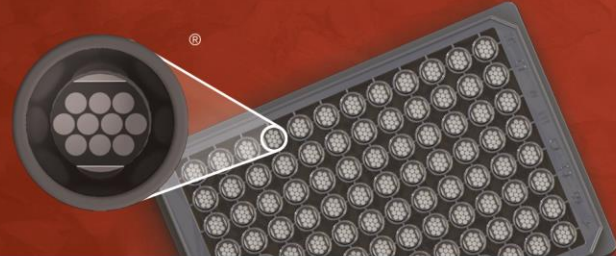


# U-PLEX<sup>®</sup> Mouse EPO Assay



[www.mesoscale.com](http://www.mesoscale.com)<sup>®</sup>

## Ordering Information

MSD<sup>®</sup> Customer Service  
 Phone: 1-301-947-2085  
 Fax: 1-301-990-2776  
 Email: [CustomerService@mesoscale.com](mailto:CustomerService@mesoscale.com)

## Scientific Support

Phone: 1-301-947-2025  
 Email: [ScientificSupport@mesoscale.com](mailto:ScientificSupport@mesoscale.com)

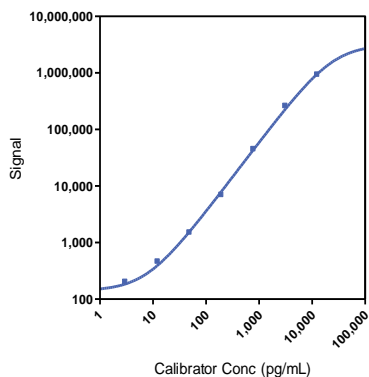
## Company Address

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

<b>Product Options</b>	Available as part of U-PLEX Biomarker Group 1 (ms) multiplex combination: K15069L-1/-2/-4
	Individual assay: K152VXK-1/-2/-4; Antibody Set: B22VX-2/B22VX-3
	For more ordering options, please visit <a href="http://www.mesoscale.com">www.mesoscale.com</a>
<b>Instrument Compatibility</b>	SECTOR <sup>®</sup> Imager 2400, SECTOR Imager 6000, MESO <sup>®</sup> SECTOR S 600, MESO QuickPlex <sup>®</sup> SQ 120
<b>Sample Type</b>	Mouse serum, EDTA plasma, and cell culture supernatants
<b>Assay Protocol</b>	Refer to the U-PLEX Biomarker Group 1 (Mouse) product insert available at <a href="http://www.mesoscale.com">www.mesoscale.com</a>

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 EPO 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<sup>™</sup> Small Spot Streptavidin plates.

## Representative Calibration Curve and Sensitivity



Assay	Median LLOD (pg/mL)	LLOD Range (pg/mL)
EPO	4.5	4.5-4.7

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 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
EPO	High	4,754	2.2	4.3
	Mid	450	2.3	5.1
	Low	50	2.8	6.8

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
EPO	High	78	72-83	54	52-59	112	102-120
	Mid	72	63-77	52	46-57	108	97-123
	Low	75	66-81	60	54-64	97	90-109

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	Spiked Serum	Spiked Plasma
Median (pg/mL)	ND	ND	218	111
Range (pg/mL)	ND-121	ND-225	218-225	110-117
% Detected	30	30	100	100

ND = non-detectable (< 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
EPO	2	120	114-129	2	99	91-105	2	93	88-101
	4	112	76-134	4	95	83-106	4	88	83-96
	8	116	75-137	8	93	80-104	8	85	78-94

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, EPO Antibody Set was tested individually against a larger panel of recombinant mouse analytes for nonspecific binding (GM-CSF, IFN- $\gamma$ , IL-1 $\beta$ , IL-2, IL-4, IL-5, IL-6, IL-10, IL-12p70, IL-13, KC/GRO, TNF- $\alpha$ , EPO, 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 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 EPO is included in Calibrator 5 blend. The full-length recombinant mouse protein expressed in murine myeloma cell line is used.

**Antibodies:** The U-PLEX Mouse EPO 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, & 10-SPOT (designs), 384 WELL 1 & 4-SPOT (designs), MSD (design), U-PLEX (design), S-PLEX (design), V-PLEX (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.

©2015 Meso Scale Diagnostics, LLC. All rights reserved.

