



[www.mesoscale.com](http://www.mesoscale.com)

## Ordering Information

MSD Customer Service  
Phone: 1-240-314-2795  
Fax: 1-301-990-2776  
Email: [CustomerService@mesoscale.com](mailto:CustomerService@mesoscale.com)

## Scientific Support

Phone: 1-240-314-2798  
Email: [ScientificSupport@mesoscale.com](mailto:ScientificSupport@mesoscale.com)

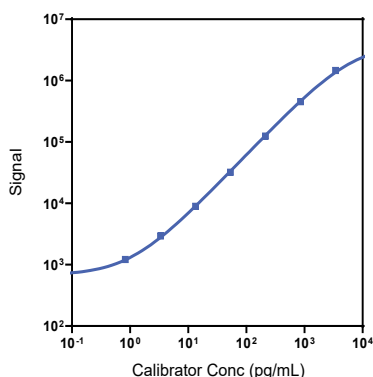
## Company Address

Meso Scale Discovery  
A division of  
Meso Scale Diagnostics, LLC.  
1601 Research Boulevard  
Rockville, MD 20850-3173 USA

Product Options	Catalog Number	Description
Multiplex	K15068M, K25068M	U-PLEX Biomarker Group 1 (NHP)
Singleplex	K156WIK-1/-2/-4	U-PLEX NHP IL-22 Assay with SECTOR™ plates
	K156WIK-21/-22/-24	U-PLEX NHP IL-22 Assay with QuickPlex Ultra™ plates
	K256WIK-2/-4	U-PLEX NHP IL-22 Assay with 384-well plates
Antibody Set	B21WI-2/-3	U-PLEX Human IL-22 Assay Antibody Set
Assay Protocol	U-PLEX Product Inserts are available at <a href="http://www.mesoscale.com">www.mesoscale.com</a>	

The MESO SCALE DISCOVERY® 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® NHP IL-22 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 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-22	0.13	0.08-0.19

The Calibrator curve was fitted with a 4-parameter logistic model with a  $1/Y^2$  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
IL-22	High	836	3.2	6.4
	Mid	98	4.4	9.6
	Low	11	4.2	11.1

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.**

## Spike Recovery

	Spike Level	Serum (N=5)		Plasma (N=5)		Cell Culture Media (N=5)	
		Average % Recovery	% Recovery Range	Average % Recovery	% Recovery Range	Average % Recovery	% Recovery Range
Cynomolgus Monkey	High	79	61-89	95	85-101	110	104-114
	Mid	72	62-79	81	68-89	114	111-119
	Low	62	52-68	72	63-77	115	112-121
Rhesus Monkey	High	81	59-94	91	83-95	110	104-114
	Mid	72	46-88	91	78-95	114	111-119
	Low	68	39-84	86	69-88	115	112-121

Normal 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.

$$\% \text{ Recovery} = (\text{measured concentration} / \text{expected concentration}) \times 100$$

## Tested Samples

	Sample Type	Serum (N=11)	Plasma (N=11)	Cell Culture Media (N=10)
Cynomolgus Monkey	Median (pg/mL)	ND	ND	3.5
	Range (pg/mL)	ND-2.0	ND-3.3	0.4-124
	% Detected	27	18	100
Rhesus Monkey	Median (pg/mL)	0.14	ND	0.8
	Range (pg/mL)	ND-428	ND-2.6	ND-14
	% Detected	54	36	80

Normal serum, EDTA plasma, and cell culture media were diluted 2-fold prior to the assay. ND = not detectable (<LLOD)

## Dilution Linearity

	Fold Dilution	Serum (N=5)		Fold Dilution	Plasma (N=5)		Fold Dilution	Cell Culture Media (N=4)	
		Average % Recovery	% Recovery Range		Average % Recovery	% Recovery Range		Average % Recovery	% Recovery Range
Cynomolgus Monkey	2	112	103-120	2	111	104-117	2	114	111-118
	4	120	104-131	4	119	108-139	4	112	108-118
	8	120	101-136	8	115	102-142	8	106	97-113
Rhesus Monkey	2	111	102-119	2	111	107-122	2	114	111-118
	4	110	99-119	4	110	100-123	4	112	108-118
	8	108	101-118	8	112	98-130	8	106	97-113

Normal 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.

$$\% \text{ Recovery} = (\text{measured concentration} / \text{expected concentration}) \times 100$$

# MSD U-PLEX NHP IL-22

## Specificity

To assess specificity, the IL-22 Antibody Set was tested individually against a larger panel of recombinant human analytes for nonspecific binding (CTACK, Eotaxin, Eotaxin-2, Eotaxin-3, ENA-78, FLT3L, Fractalkine, G-CSF, GM-CSF, GRO- $\alpha$ , I-309, IFN- $\alpha$ 2a, IFN- $\gamma$ , IL-1 $\alpha$ , IL-1 $\beta$ , IL-1RA, IL-2, IL-4, IL-5, IL-6, IL-7, IL-8, IL-9, IL-10, IL-12/IL-23p40, IL-12p70, IL-13, IL-15, IL-16, IL-17A, IL-17AF, IL-17B, IL-17C, IL-17D, IL-17F, IL-18, IL-22, IL-23, IP-10, I-TAC, MCP-1, MCP-2, MCP-3, MCP-4, M-CSF, MDC, MIF, MIP-1 $\alpha$ , MIP-1 $\beta$ , MIP-3 $\alpha$ , MIP-3 $\beta$ , MIP-5, SDF-1 $\alpha$ , TARC, TNF- $\alpha$ , TNF- $\beta$ , TPO, TRAIL, VEGF-A, and YKL-40). Nonspecific binding was less than 0.5%.

$$\% \text{ Nonspecificity} = (\text{nonspecific signal} / \text{specific signal}) \times 100$$

## 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 your assay needs, other diluents may be tested.

## Assay Components

**Calibrator:** IL-22 is included in Calibrator 6. The full-length recombinant protein is expressed in *E. coli*.

**Antibodies:** The U-PLEX NHP IL-22 Assay uses a rat monoclonal antibody for capture and a 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.

