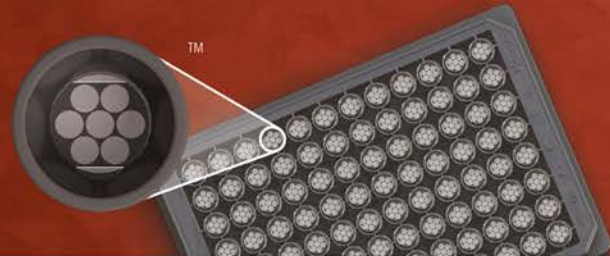


MSD[®] Rat Demonstration 7-Plex Ultra-Sensitive Kit

For quantitative determination in rat serum and plasma



Alzheimer's Disease
BioProcess
Cardiac
Cell Signaling
Clinical Immunology
Cytokines
Hypoxia
Immunogenicity
Inflammation
Metabolic
Oncology
Toxicology
Vascular

Catalog Numbers

Rat Demonstration 7-Plex Ultra-Sensitive Kit	
Kit size	
1 plate	K15014C-1
5 plates	K15014C-2
25 plates	K15014C-4

Ordering information

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

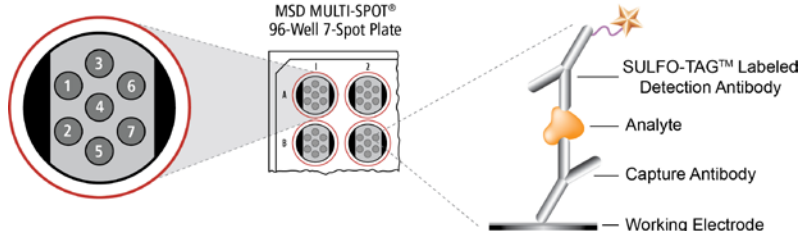
Company Address

MESO SCALE DISCOVERY[®]
A division of
Meso Scale Diagnostics, LLC.
9238 Gaither Road
Gaithersburg, MD 20877 USA

www.mesoscale.com[®]

For Research Use Only.
Not for use in diagnostic procedures.

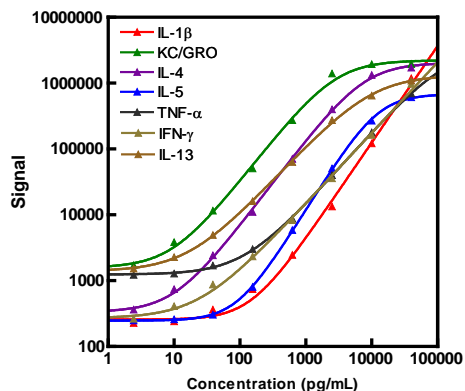
- | | |
|------------------|-------------------|
| 1 = IL-1 β | 5 = TNF- α |
| 2 = KC/GRO/CINC | 6 = IFN- γ |
| 3 = IL-4 | 7 = IL-13 |
| 4 = IL-5 | |



Cytokines are immunomodulating agents mediating a number of physiological responses, including immunity, inflammation, and a variety of pathophysiological conditions. In disease states, the circulating levels of cytokines are altered, making them valuable biomarkers that can be associated with many clinical conditions. MESO SCALE DISCOVERY[®] offers a wide range of cytokine assays in singleplex and multiplex formats. The Rat Demonstration 7-Plex Ultra-Sensitive Kit can be used to simultaneously assess the levels of 7 cytokines in serum and plasma samples; it may also be used with bronchial lavage, wound fluids, urine, and other biological samples. Tissue culture kits are available for cell supernatant samples. Visit www.mesoscale.com for a complete list of human, rodent, canine, bovine, and non-human primate assays.

Assay Sensitivity

The following standard curves illustrate the dynamic range of the assays in the Rat Demonstration 7-Plex Ultra-Sensitive Kit.



	IL-1 β	KC/GRO/CINC	IL-4	IL-5
LLOD (pg/mL)	25	1.3	1.6	25
	TNF- α	IFN- γ	IL-13	
LLOD (pg/mL)	23	9.0	2.3	

The lower limit of detection (LLOD) is a calculated concentration based on a signal 2.5 standard deviations above the blank (zero) calibrator. The values shown represent the average LLOD over multiple kit lots.

MSD Advantages

- **Multiplexing:** Multiple analytes can be measured in one well using typical sample volumes of 25 μ L or less without compromising speed or performance
- **Large dynamic range:** Linear range of up to five logs enables the measurement of native levels of biomarkers in normal and diseased samples without multiple dilutions
- **Minimal background:** The stimulation mechanism (electricity) is decoupled from the signal (light)
- **Simple protocols:** Only labels near the electrode surface are detected, enabling assays with fewer washes
- **Flexibility:** Labels are stable, non-radioactive, and conveniently conjugated to biological molecules
- **High sensitivity and precision:** Multiple excitation cycles of each label enhance light levels and improve sensitivity

For a complete list of products, please visit our website at www.mesoscale.com

MSD Cytokine Assays

Specificity

Cross-reactivity of the assays within the same well was measured using a single analyte at 10 000 or 40 000 pg/mL. The high values were chosen to yield a high specific signal and to provide a stringent test for cross-reactivity. The signal measured for the other analytes were used to calculate cross-reactivity. The % cross-reactivity was less than 0.5% for all analytes.

	% Specificity						
	IL-1 β	KC/GRO	IL-4	IL-5	TNF- α	IFN- γ	IL-13
IL-1 β	100	0.12	0.12	0.09	0.04	0.08	0.02
KC/GRO	0.01	100	<0.01	<0.01	<0.01	<0.01	<0.01
IL-4	0.20	0.08	100	0.15	0.08	0.19	0.04
IL-5	0.08	0.11	0.08	100	0.06	0.12	0.04
TNF- α	0.42	0.17	0.24	0.37	100	0.11	0.35
IFN- γ	0.08	0.06	0.10	0.15	0.24	100	0.06
IL-13	<0.01	<0.01	<0.01	<0.01	<0.01	0.03	100

Precision

Control samples were spiked with high, mid, and low levels of each analyte and were run in triplicate over multiple days ($n>3$) using multiple plate lots. Average Intra-plate %CV is the average %CV of the control replicates within an individual plate. Inter-plate %CV expresses the variability of controls across 13 plates over multiple days.

	Level	%CV						
		IL-1 β	KC/GRO	IL-4	IL-5	TNF- α	IFN- γ	IL-13
Average Intra-Plate	Low	5.3	2.2	2.6	5.6	3.3	3.3	3.0
	Mid	4.5	2.7	5.4	2.3	3.7	3.5	2.6
	High	4.2	6.0	2.5	4.3	3.7	5.9	4.9
Inter-Plate	Low	8.9	9.2	6.1	9.5	8.0	9.6	10.5
	Mid	8.7	8.4	7.8	9.1	6.0	7.7	8.6
	High	5.4	7.9	9.3	5.4	9.9	7.1	7.1

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, WWW.MESOSCALE.COM, MSD, MSD (DESIGN), DISCOVERY WORKBENCH, QUICKPLEX, MULTI-ARRAY, MULTI-SPOT, SULFO-TAG, SECTOR, SECTOR HTS, SECTOR PR, 7-SPOT (DESIGN) and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.
© 2012 Meso Scale Diagnostics, LLC. All rights reserved.

For Research Use Only. Not for use in diagnostic procedures.

