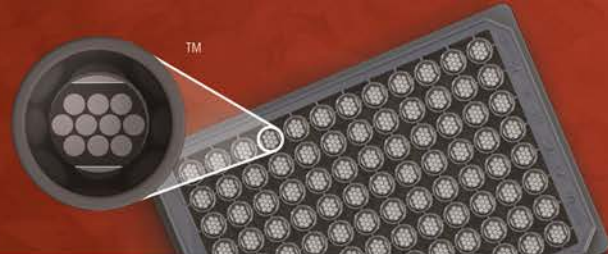


MSD® Mouse TH1/TH2 9-Plex Ultra-Sensitive Kit

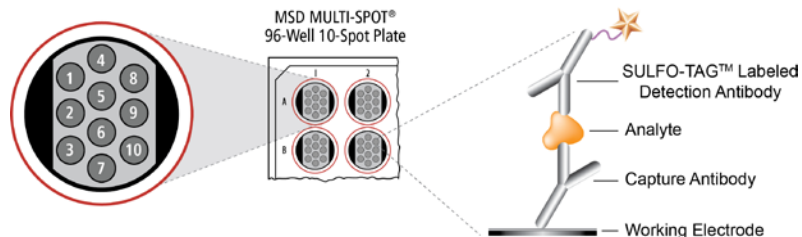
For quantitative determination in mouse serum and plasma



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

Catalog Numbers

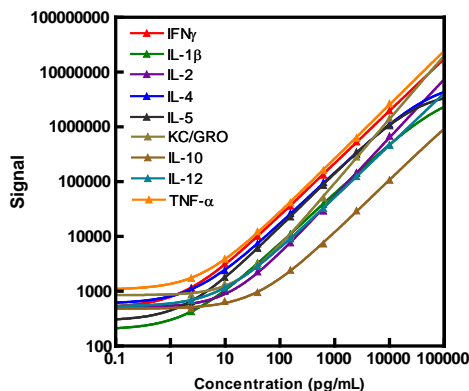
- | | |
|-------------------|------------------------|
| 1 = IFN- γ | 6 = KC/GRO/CINC |
| 2 = IL-1 β | 7 = IL-10 |
| 3 = IL-2 | 8 = IL-12 total |
| 4 = IL-4 | 9 = <i>BSA Blocked</i> |
| 5 = IL-5 | 10 = TNF- α |



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 Mouse TH1/TH2 9-Plex Ultra-Sensitive Kit can be used to simultaneously assess the levels of 9 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 Mouse TH1/TH2 9-Plex Ultra-Sensitive Kit.



	IFN- γ	IL-1 β	IL-2	IL-4	IL-5
LLOD (pg/mL)	0.47	2.1	3.0	0.87	0.70

	KC/GRO/CINC	IL-10	IL-12 total	TNF- α
LLOD (pg/mL)	2.9	11	5.3	1.0

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.

Ordering information

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

Mouse TH1/TH2 9-Plex Ultra-Sensitive Kit	
Kit size	
1 plate	K15013C-1
5 plates	K15013C-2
25 plates	K15013C-4

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

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, 10-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.