



MSD[®] Bioanalytical Laboratory Validated Assay List

V-PLEX® Testing Service — Proinflammatory Panel 1 Human (S005CD-1)

Analyte	LLOD (pg/mL)	LLOQ (pg/mL)	ULOQ (pg/mL)	Dynamic Range (pg/mL)
IFN-γ	0.37	1.76	938	0.37 – 938
IL-1β	0.05	0.646	375	0.05 – 375
IL-2	0.09	0.89	938	0.09 – 938
IL-4	0.02	0.218	158	0.02 – 158
IL-6	0.06	0.633	488	0.06 – 488
IL-8	0.07	0.591	375	0.07 – 375
IL-10	0.04	0.298	233	0.04 – 233
IL-12p70	0.11	1.22	315	0.11 – 315
IL-13	0.24	4.21	353	0.24 – 353
TNF-α	0.04	0.69	248	0.04 – 248

Validated* in serum and EDTA plasma

V-PLEX Testing Service — Cytokine Panel 1 Human (S007CD-1)

Analyte	LLOD (pg/mL)	LLOQ (pg/mL)	ULOQ (pg/mL)	Dynamic Range (pg/mL)
GM-CSF	0.16	0.842	750	0.16 – 750
IL-1α	0.09	2.85	278	0.09 – 278
IL-5	0.14	4.41	562	0.14 – 562
IL-7	0.12	0.851	563	0.12 – 563
IL-12/IL-23p40	0.33	1.32	2,250	0.33 – 2,250
IL-15	0.15	0.774	525	0.15 – 525
IL-16	2.83	19.1	1,870	2.83 – 1,870
IL-17A	0.31	3.19	3,650	0.31 – 3,650
TNF-β	0.08	0.465	458	0.08 – 458
VEGF-A	1.12	7.70	562	1.12 – 562

Validated* in serum and EDTA plasma

V-PLEX Testing Service — Chemokine Panel 1 Human (S006CD-1)

Analyte	LLOD (pg/mL)	LLOQ (pg/mL)	ULOQ (pg/mL)	Dynamic Range (pg/mL)
Eotaxin	3.26	12.3	1,120	3.26 – 1,120
Eotaxin-3	1.77	10.2	3,750	1.77 – 3,750
IL-8 (HA)	95.6	713	43,400	95.6 – 43,400
IP-10	0.37	1.37	500	0.37 – 500
MCP-1	0.09	1.09	375	0.09 – 375
MCP-4	0.18	1.49	472	0.18 – 472
MDC	1.22	88.3	3,700	1.22 – 3,700
MIP-1α	3.02	13.8	743	3.02 – 743
MIP-1β	0.17	1.88	520	0.17 – 520
TARC	0.22	3.32	1,120	0.22 – 1,120

Validated* in serum and EDTA plasma

Custom Sample Testing — U-PLEX® Interferon Combo Human (S01CD-1)

Analyte	LLOD (pg/mL)	LLOQ (pg/mL)	ULOQ (pg/mL)	Dynamic Range (pg/mL)
IFN-γ	7.4	32	29,250	7.42 – 29,250
IFN-α2a	1.8	8.8	32,400	1.77 – 32,400
IFN-β	7.8	53	97,200	7.77 – 97,200
IL-29/IFN-λ1	1.4	4.2	7,686	1.38 – 7,685

Validated* in serum and EDTA plasma

S-PLEX® Testing Service

Analyte	LLOD (fg/mL)	LLOQ (fg/mL)	ULOQ (fg/mL)	Dynamic Range (fg/mL)
Human Tau (pT217) (S035CD-1)	395	1,840	3,760,000	395 – 3,760,000
Human Tau (pT181) (S036CD-1)	84.7	1,080	2,210,000	84.7 – 2,210,000

Validated* in serum, EDTA plasma, heparin plasma, and CSF

S-PLEX Testing Service — Proinflammatory Panel 1 Human (S038CD-1)

Analyte	LLOD (fg/mL)	LLOQ (fg/mL)	ULOQ (fg/mL)	Dynamic Range (fg/mL)
IFN-γ	9.0	127	25,400	9.0 – 25,400
IL-10	19	174	139,000	19 – 139,000
IL-12p70	43	972	389,000	43 – 389,000
IL-1β	71	672	134,000	71 – 134,000
IL-2	12	321	64,200	12 – 64,200
IL-4	8.0	54	43,400	8.0 – 43,400
IL-6	11	108	43,120	11 – 43,120
TNF-α	11	105	42,100	11 – 42,100
IL-17A	44	2,070	207,000	44 – 207,000

Validated* in serum and EDTA plasma

*Analytical method validation includes accuracy, precision, sensitivity, dilution linearity, and selectivity testing performed over multiple testing days by multiple operators.

Custom Assays and Services

- Custom panels can be developed and validated as U-PLEX assays
- Additional qualification or validation — from exploratory to full method validation — available upon request
- Analytes in any panel may be selected and analyzed as a singleplex
- Qualifications may be customized based on the context of use
- Matrix testing beyond currently validated matrices also available
- Contact MSD Customer Service for more information about developing and validating custom assays

Highlights

- Expert MSD Bioanalytical Lab analysts who can support GLP and GCLP studies
- Established Quality Management System that complies with the applicable principles of GLP and GCLP
- Assays may be qualified for basic research or designed for a full method validation
- Fit-for-purpose custom validated assays to support a comprehensive bioanalytical strategy including:
 - Pharmacokinetic assays
 - Screening anti-drug (ADA) assays
 - Confirmatory ADA assays
 - Neutralizing ADA assays (cell-based and ligand binding assays)
 - Serology Assays
- Seamless analytical method transfer after method development for expedited validation and/or sample analysis
- Reduced variability due to specific kit builds encompassing validation and sample analysis, eliminating need for lot to lot qualification
- Assays developed in standard (e.g., V-PLEX, S-PLEX) and novel (e.g., N-PLEX®) formats that are not readily available via CROs

Frequently Asked Questions

Q. When will I receive the data?

A. Our goal is to provide your data within five business days of receipt of your samples.

Q. How will I get the data?

A. You will receive the data via email to a designated contact or recipient. The data file can be imported into your LIMS or other data analysis systems.

Q. What if I anticipate having two or more batches of samples?

A. There are two options: a) we can wait until all batches of samples are received to run the most efficient plate; or b) we can run each batch of samples as we receive them. In this second option, you will need to purchase additional plates.

Q. How are samples run?

A. We can run up to 30 samples in duplicate on each plate. For more than 30 samples, you can request additional plates. If you have fewer than 30 samples, you can request more replicates per sample or leave the empty wells blank.

Q. How much sample volume do I need?

A. You will need a minimum of 100 µL of sample for one measurement in duplicate.

Q. Can you perform a dilution series?

A. Yes. Each point in the dilution series is counted as an individual sample.

Q. Can you handle special matrices and/or sample processing steps?

A. Yes. We are happy to discuss your sample types and processing requirements to understand your needs. Depending on the complexity, certain processes can be done at no additional charge, and others may incur costs. Our team has experience with varied sample types and matrices.

Q. Are there different requirements for the samples for running V-PLEX or S-PLEX assays?

A. No.

Q. How do I send my samples?

A. Please contact us at Bioanalytical.Services@mesoscale.com. We will send you a form to accompany your shipment. The shipping address is included on the form.

Ask us how we can help;
Schedule a call with our team:

Phone: 1-240-314-2795

Fax: 1-301-990-2776

Email: CustomerService@mesoscale.com

Bioanalytical.Services@mesoscale.com

MESO SCALE DISCOVERY, Meso Scale Diagnostics, MSD, mesoscale.com, www.mesoscale.com, methodicalmind.com, www.methodicalmind.com, DISCOVERY WORKBENCH, InstrumentLink, MESO, MesoSphere, Methodical Mind, MSD GOLD, MULTI-ARRAY, MULTI-SPOT, QuickPlex, ProductLink, SECTOR, SECTOR HTS, SECTOR PR, SULFO-TAG, TeamLink, TrueSensitivity, TURBO-BOOST, TURBO-TAG, N-PLEX, R-PLEX, S-PLEX, T-PLEX, U-PLEX, V-PLEX, MSD (design), MSD (luminous design), Methodical Mind (head logo), 96 WELL SMALL-SPOT (design), 96 WELL 1-, 4-, 7-, 9-, & 10-SPOT (designs), 384 WELL 1- & 4-SPOT (designs), N-PLEX (design), R-PLEX (design), S-PLEX (design), T-PLEX (design), U-PLEX (design), V-PLEX (design), It's All About U, Spot the Difference, The Biomarker Company, and The Methodical Mind Experience are trademarks and/or service marks owned by or licensed to Meso Scale Diagnostics, LLC. All other trademarks and service marks are the property of their respective owners.

© 2023 Meso Scale Diagnostics, LLC. All rights reserved.



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