



MSD[®] Biomarker Services

Whether you require a biomarker screen or the development of a validated kit for lead biomarkers, MESO SCALE DISCOVERY[®] Biomarker Services will tailor a custom service package for you. With over twenty years of experience in developing and multiplexing high-quality commercial immunoassays, MSD can accelerate your biomarker program from early stage discovery to the production of fully validated kits or components.

SCREEN

CRP	IL-6	MDC
Eotaxin	IL-7	MIP-1 α
Eotaxin-3	IL-8	MIP-1 β
FGF (basic)	IL-10	PIGF
Flt-1	IL-12/IL-23p40	SAA
GM-CSF	IL-12p70	TARC
ICAM-1	IL-13	Tie-2
IFN- γ	IL-15	TNF- α
IL-1 α	IL-16	TNF- β
IL-1 β	IL-17	VCAM-1
IL-2	IP-10	VEGF-A
IL-4	MCP-1	VEGF-C
IL-5	MCP-4	VEGF-D

MSD screens your samples in multiplex to identify potential biomarkers

Over 135 high-impact analytes measured with a rapid turnaround time

Great for limited sample volumes

SELECT

<input type="checkbox"/> CRP	<input type="checkbox"/> IL-1 β
<input type="checkbox"/> Eotaxin	<input type="checkbox"/> IL-2
<input checked="" type="checkbox"/> Eotaxin-3	<input checked="" type="checkbox"/> IL-4
<input type="checkbox"/> FGF (basic)	<input type="checkbox"/> IL-5
<input type="checkbox"/> Flt-1	<input type="checkbox"/> IL-6
<input checked="" type="checkbox"/> GM-CSF	<input checked="" type="checkbox"/> IL-7
<input type="checkbox"/> ICAM-1	<input type="checkbox"/> IL-8
<input checked="" type="checkbox"/> IFN-γ	<input type="checkbox"/> IL-10
<input type="checkbox"/> IL-1 α	<input type="checkbox"/> IL-12/IL-23p40

MSD develops customized assays for your lead biomarkers from a screen

Fit-for-purpose assays designed for optimal test development

Transition from identification of promising biomarkers to final panels

VALIDATE

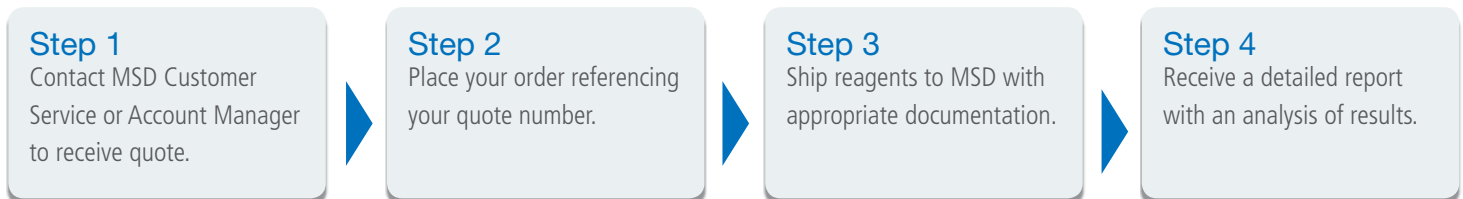


MSD develops and validates your final biomarker panel

Manufactured under ISO 9001:2008 certified quality management system

Your validated biomarker test is available for order

Contact us to discuss your sample testing service needs.



Phone: 1-240-314-2793 . Fax: 1-301-990-2776
Email: assayservices@mesoscale.com

Human Biomarker Screening Service 1 (Catalog # S01CA-1/-2/-3)

A1AT NEW	E-Cadherin	IFN- α	IL-17E/IL-25 NEW	MIG	Resistin
A2M	ENA-78	IFN- γ	IL-17F NEW	MIP-1 α	SAA
Active GLP-1	Endoglin	IL-1 α	IL-18	MIP-1 β	SCF
Adiponectin	Eotaxin	IL-1 β	IL-21	MIP-3 α	SDF-1 α
Angiopoietin 1	Eotaxin-2	IL-1RA	IL-22 NEW	MIP-3 β	TARC
Angiopoietin 2	Eotaxin-3	IL-2	IL-23 NEW	MIP-5	Tenascin C
APOC3 NEW	EPO	IL-4	IL-27 NEW	MMP-1	Thrombomodulin
B2M	E-Selectin	IL-5	IL-29/IFN- λ 1 NEW	MMP-3	Tie-2
Calbindin	FABP3	IL-6	IL-31 NEW	MMP-9	TIMP-1 NEW
CA 15.3	Factor VII	IL-6R	IL-33	Myl3	TNF- α
CA 50	Fas	IL-7	Insulin	Myoglobin	TNF- β
CA 125	FasL	IL-8	IP-10	Nectin-4	TNF-RI
CEA	FGF (basic)	IL-10	I-TAC	NT-proBNP	TNF-RII
c-Kit	Flt-1	IL-12/IL-23p40	Leptin	Osteoactivin	TPO
CKMB	Flt-3 Ligand	IL-12p70	MCP-1	Osteocalcin	TRAIL
Clusterin	Fractalkine	IL-13	MCP-2	Osteonectin	TSLP NEW
C-Peptide	GIP	IL-15	MCP-3	Osteopontin	VCAM-1
CRP	Glucagon	IL-16	MCP-4	Osteoprotegerin	VEGF-A
CTACK	GM-CSF	IL-17A	M-CSF	P-Cadherin	VEGF-C
cTnI	GRO- α	IL-17A (Gen B) NEW	MDC	PIGF	VEGF-D
cTnT	I-309	IL-17A/F NEW	Mesothelin	P-Selectin	YKL-40
Cystatin C NEW	ICAM-1	IL-17B	Met	PYY (total)	
Cytokeratin-8	ICAM-3	IL-17D	MIF	RANTES	

Human Biomarker Screening Service 2 (Catalog # S02CA-1/-2/-3) **NEW**

Eotaxin	IL-4	IL-13	IL17E/IL-25	IL-31	OPGN
Eotaxin-3	IL-5	IL-15	IL-17F	IL-33	TARC
GM-CSF	IL-6	IL-16	IL-18	IP-10	TNF- α
IFN- α	IL-7	IL-17A	IL-21	MCP-1	TNF- β
IFN- γ	IL-8	IL-17A (Gen B)	IL-22	MCP-4	TSLP
IL-1 α	IL-10	IL-17A/F	IL-23	MDC	VEGF-A
IL-1 β	IL-12/IL-23p40	IL-17B	IL-27	MIP-1 α	
IL-2	IL12p70	IL-17D	IL-29/IFN- λ 1	MIP-1 β	

Human V-PLEX[®] Screening Service 1 (Catalog # S01CB-1/-2/-3)

CRP	IFN- γ	IL-7	IL-16	MIP-1 β	VCAM-1
Eotaxin	IL-1 α	IL-8	IL-17A	PIGF	VEGF-A
Eotaxin-3	IL-1 β	IL-10	IP-10	SAA	VEGF-C
FGF (basic)	IL-2	IL-12/IL-23p40	MCP-1	TARC	VEGF-D
Flt-1	IL-4	IL-12p70	MCP-4	Tie-2	
GM-CSF	IL-5	IL-13	MDC	TNF- α	
ICAM-1	IL-6	IL-15	MIP-1 α	TNF- β	

Results will be provided in Microsoft Excel format and include assay range, limit of detection, and calculated concentrations derived for each of the samples tested.

MESO SCALE DISCOVERY, MSD, V-PLEX, www.mesoscale.com, SPOT THE DIFFERENCE, and MSD (design) are trademarks and/or service marks of Meso Scale Diagnostics, LLC. © 2016 Meso Scale Diagnostics, LLC. All rights reserved.



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

MK-D5-221-v2-2016/June