

# R-PLEX<sup>®</sup> Antibody Set Dilution Factors and Assay/Antibody Diluent Combinations

The tables below provide the dilution factors and primary diluent combinations for samples tested with R-PLEX Antibody Sets. These dilution factors and diluent combinations are recommended as starting points when developing your assay. However, other sample dilutions and diluents should be considered based on the sample environment. Please refer to the R-PLEX datasheet for additional information on your analytes of interest. For further guidance, please contact our Scientific Support team at [scientificsupport@mesoscale.com](mailto:scientificsupport@mesoscale.com).

## Notes:

- (1) Additional assay diluent is necessary for samples that are diluted greater than 10-fold. Diluent 100 may be used in place of the assay diluent for samples that require high dilution. Testing of different diluents can help optimize assays for specific experimental conditions.
- (2) Some assay and antibody diluents require additives, as noted in the tables. For more information, refer to the product-specific datasheet available at [www.mesoscale.com/R-PLEX-documents](http://www.mesoscale.com/R-PLEX-documents).
- (3) For cell lysate samples, the fold dilution or total protein amount ( $\mu\text{g}/\text{well}$ ) is dependent on the cell type and model system used. We recommend performing a dilution series and testing for diluent compatibility.

Human							
Antibody Set	Dilution Factor					Primary Assay Diluent	Primary Antibody Diluent
	Serum	Plasma	Urine	CSF	Other		
4-1BB/TNFRSF9	Neat	Neat	—	—	—	12	11
4-1BBL/TNFSF9	Neat	Neat	—	—	—	12	11
5'-Nucleotidase	2	2	—	—	—	12	11
6Ckine/CCL21	2	2	—	—	—	57	3
A2M	4,000	4,000	—	—	—	12	12
A $\beta$ (total)	—	—	—	4	—	35	100
A $\beta$ 38 (4G8)	—	—	—	2	—	35	100
A $\beta$ 38 (6E10)	—	—	—	2	—	35	100
A $\beta$ 40 (4G8)	—	—	—	2	—	35	100
A $\beta$ 40 (6E10)	—	—	—	3	—	35	100
A $\beta$ 42 (4G8)	—	—	—	2	—	35	100
A $\beta$ 42 (6E10)	—	—	—	2	—	35	100
ACE	2	2	—	—	—	57	3
ACE2	Neat	Neat	—	—	—	12	12
AChE	Neat	Neat	—	—	—	57	3
ADAM 9	Neat	Neat	—	—	—	12	11
ADAM17	Neat	Neat	—	—	—	12	11
ADAM-TS 13	4	4	—	—	—	57	3
ADP-ribosyl cyclase	Neat	Neat	—	—	—	57	3
AFP	Neat	Neat	—	—	—	57	3
AGP	—	300,000	—	—	—	37	37
$\alpha$ -GST	2	2	—	—	—	37	37
Aiolos	—	—	—	—	0.62 $\mu\text{g}/\text{well}$ (cell lysate)	39 (with additives)	MSD Tris Wash Buffer (with additives)



## R-PLEX Antibody Set Dilution Factors and Assay/Antibody Diluent Combinations

Human							
Antibody Set	Dilution Factor					Primary Assay Diluent	Primary Antibody Diluent
	Serum	Plasma	Urine	CSF	Other		
Albumin	320,000	320,000	500	—	—	100	59
ALCAM/CD166	64	64	—	—	—	57	3
Alpha-amylase 1	Neat	Neat	—	—	—	58	3
Alpha Defensin 1	8	8	—	—	—	57	3
AMICA1	2	2	—	—	—	13	11
Amphiregulin	Neat	Neat	—	—	—	57	3
Ang-1	2	2	2	2	—	10	3
Ang-2	2	2	2	2	—	10	3
Ang-4	Neat	Neat	—	—	—	57	3
Angiogenin	10,000	10,000	—	—	—	57	3
Ang-like 3	25	25	—	—	—	13	11
Ang-like 4	Neat	Neat	—	—	—	13	11
Ang-like 6	Neat	Neat	—	—	—	13	11
Annexin A1	—	4	—	—	—	7	37
Annexin A5	Neat	Neat	—	—	—	13	11
Antileukoproteinase	50	50	—	—	—	35	35
ApoA1	—	300,000	—	—	—	37	37
ApoC3	50,000	50,000	—	—	—	101	100
ApoE	100	100	—	—	—	57	3
APRIL/TNFSF13	8	—	—	—	—	57	3
Arginase-1	2	2	—	—	—	58	3
Autotaxin	16	8	—	—	—	57	3
AXL/UFO	8	8	—	—	—	57	3
B2M	4,000	4,000	4,000	4,000	—	12	12
B7-H6	Neat	Neat	—	—	—	58	3
Basigin	4	4	—	—	—	12	11
BCA-1/BLC	2	2	—	—	—	57	3
Bcl-2	—	—	—	—	0.31 µg/well (cell lysate)	39 (with additives)	39
Bcl-X/Bcl2-L-1	—	—	—	—	0.31 µg/well (cell lysate)	39 (with additives)	39
Bikunin	4,000	4,000	—	—	—	57	3
BIM/Bcl2-L-11	—	—	—	—	2.0 µg/well (cell lysate)	39 (with additives)	39
BMP-2	Neat	Neat	—	—	—	7	11
BMP-4	Neat	Neat	—	—	—	7	11
BMP-5	Neat	Neat	—	—	—	57	3
BMP-6	Neat	Neat	—	—	—	57	3
BMP-7	Neat	Neat	—	—	—	7	11
BMP-9/GDF-2	8	8	—	—	—	7	11
BMP-11/GDF-11	Neat	Neat	—	—	—	12	11
C1q receptor	1,000	1,000	—	—	—	57	3
CA1	—	5,000	—	—	—	101	37
CA15-3	20	20	20	—	—	58	3
CA125	20	20	20	20	—	58	3
Calbindin	10	10	10	—	10	7	3
Calprotectin	8	—	—	—	—	101	37
Cathepsin B (total)	4	4	—	—	—	12	11
Cathepsin D	10	10	—	—	—	101	37



Dash (—) indicates that the information for that sample type is not available.

18199-v15-2023Jun | 2

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES

## R-PLEX Antibody Set Dilution Factors and Assay/Antibody Diluent Combinations

Human							
Antibody Set	Dilution Factor					Primary Assay Diluent	Primary Antibody Diluent
	Serum	Plasma	Urine	CSF	Other		
Cathepsin L2	Neat	Neat	—	—	—	12	11
Cathepsin S (total)	16	16	—	—	—	12	11
CCN1	Neat	Neat	—	—	—	57	3
CD5	—	Neat	—	—	—	58	3
CD5L	5,000	5,000	8	—	—	58	3
CD8α	8	8	—	—	—	57	3
CD9 (EV)	10	50	4	10	—	53	53
CD14	—	—	—	—	200 (synovial fluid)	101	3
CD21/CR2	500	500	—	—	—	57	3
CD22	Neat	Neat	—	—	—	57	3
CD30/TNFRSF8	Neat	Neat	—	—	—	57	3
CD31/PECAM-1	20	20	—	—	—	12	12
CD40/TNFRSF5	2	2	—	—	—	58	3
CD44	Neat	Neat	—	—	—	12	11
CD47	Neat	Neat	—	—	—	57	3
CD63 (EV)	10	20	Neat	10	—	53	53
CD80/B7-1	Neat	Neat	—	—	—	7	11
CD81 (EV)	4	10	4	10	—	53	53
CD163	2	2	—	—	—	57	3
CD177	2	4	—	—	—	57	3
CD200	Neat	Neat	—	—	—	57	3
CEACAM-1	4	4	—	—	—	12	11
CEACAM-5	—	4	—	—	—	7	37
CEACAM-7/CGM2	Neat	Neat	—	—	—	13	11
Chemerin	100	100	—	—	—	57	3
Cholinesterase	1000	1000	—	—	—	57	3
Chromogranin-A	Neat	Neat	—	—	—	57	3
cIAP1	—	—	—	—	2.5 µg/well (cell lysate)	MSD Tris Lysis Buffer (with additives)	MSD Tris Wash Buffer
cIAP2	—	—	—	—	2.5 µg/well (cell lysate)	MSD Tris Lysis Buffer (with additives)	MSD Tris Wash Buffer
CKBB	2	2	—	—	—	101	3
CK Beta 8-1/CCL23	Neat	Neat	—	—	—	2	3
Clusterin	—	5,000	—	—	—	101	37
CNTF	Neat	Neat	—	—	—	12	11
COMP/TSP-5	32	32	—	—	—	7	11
Complement C3	—	300,000	—	—	—	37	37
Complement C3a	10,000	10,000	16	—	—	58	3
Complement C5a	Neat	Neat	—	—	—	13	11
Complement C9	—	5,000	—	—	—	101	37
Complement factor D	1,000	1,000	—	—	—	13	12
Contactin-2/TAG-1	2	2	—	—	—	13	11
Corin/ATC2	2	2	—	—	—	12	12
COX-2/PTGS2	Neat	Neat	—	—	—	57	3
CRP	1,000	1,000	5	—	—	101	101
Cystatin B	Neat	Neat	—	—	—	13	11
Cystatin C	2,000	2,000	—	—	—	12	12



Dash (—) indicates that the information for that sample type is not available.

## R-PLEX Antibody Set Dilution Factors and Assay/Antibody Diluent Combinations

Human							
Antibody Set	Dilution Factor					Primary Assay Diluent	Primary Antibody Diluent
	Serum	Plasma	Urine	CSF	Other		
Cytokeratin-8	2	2	2	2	—	10	3
Cytokeratin-19	Neat	Neat	—	—	—	57	3
Decorin	16	16	—	—	—	12	11
Delta-like 1	2	2	—	—	—	13	11
DKK-1	2	2	—	—	—	MSD Tris Wash Buffer	MSD Tris Wash Buffer
DKK-3	50	50	—	—	—	57	3
DKK-4	2	2	—	—	—	13	11
DPPIV	—	5,000	—	—	—	101	37
E-Cadherin	20	20	20	20	—	58	3
EGF	10	—	500	—	—	37	37
EGFR	64	64	—	—	—	35	35
Elastase-2	2	2	—	—	—	12	11
Endoglin	50	50	50	50	—	57	3
Endostatin	50	50	—	—	—	57	3
Endothelin-1	Neat	Neat	—	—	—	57	3
Enolase 2	5	5	—	5	—	7	7
eNOS	Neat	Neat	—	—	—	12	11
Ep-CAM	Neat	Neat	—	—	—	12	11
ER- $\alpha$	Neat	Neat	—	—	20 $\mu$ g/well (cel lysate)	39	39
ErbB2	8	8	—	—	—	12	11
ErbB3/HER3	4	4	—	—	—	12	11
ERK1/2 (total)	Neat	Neat	—	—	20 $\mu$ g/well (cell lysate)	39	39
E-Selectin	2	2	2	2	—	10	3
FABP2/I-FABP	Neat	Neat	—	—	—	57	3
FABP3/H-FABP	2	2	—	—	—	12	3
FABP4/A-FABP	2	2	—	—	—	13	11
Factor VII	100	100	—	—	—	12	12
FAP- $\alpha$ /SEPR	—	100	—	—	—	101	37
Fas (soluble)	50	50	50	50	—	7	8
FasL	2	2	2	2	—	10	3
Fc-epsilon RII	2	2	—	—	—	13	11
FGF (acidic)	Neat	Neat	—	—	—	7	11
Ferritin	500	500	—	—	1 ng/mL (tissue homogenate)	12	12
FGF (basic)	2	2	2	2	—	7	11
FGF-7	Neat	Neat	—	—	—	13	11
FGF-19	2	2	—	—	—	57	3
FGF-BP1	Neat	Neat	—	—	—	12	11
Fibronectin	100,000	100,000	—	—	—	13	11
Folate R alpha	Neat	Neat	—	—	—	57	3
Follistatin	Neat	Neat	—	—	—	57	3
GAD2/GAD-65	Neat	Neat	—	—	—	12	11
Galectin-1	Neat	Neat	—	—	—	12	11
Galectin-3	50	50	—	—	—	58	3
GAS-6/AXL-L	25	25	—	—	—	13	11
GCPII/PSMA	Neat	Neat	—	—	—	12	11



Dash (—) indicates that the information for that sample type is not available.

18199-v15-2023Jun | 4

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES

## R-PLEX Antibody Set Dilution Factors and Assay/Antibody Diluent Combinations

Human							
Antibody Set	Dilution Factor					Primary Assay Diluent	Primary Antibody Diluent
	Serum	Plasma	Urine	CSF	Other		
GDF-8	Neat	Neat	—	—	—	12	11
GDF-15	—	100	—	—	—	101	37
Gelsolin	—	100	—	—	—	101	37
GFAP	—	2	—	—	—	101	27
Glicentin	—	2 (P800)	—	—	—	12	11
Granulysin	2	2	—	—	—	12	11
GRO-β/CXCL2	Neat	Neat	—	—	—	12	12
Growth Hormone	Neat	Neat	—	—	—	57	3
GRP	Neat	Neat	—	—	—	12	11
Haptoglobin	—	300,000	—	—	—	37	37
HAVCR1/KIM-1	—	—	10	—	—	37	37
HB-EGF	2	2	—	—	—	57	3
HCC-4/CCCL16	50	50	—	—	—	13	11
HE4/WFDC2	4	4	—	—	—	13	11
Heme Oxygenase 1	Neat	Neat	—	—	—	13	11
Hemoglobin alpha	400	400	—	—	10 (stool)	100	100
Hemojuvelin	Neat	Neat	—	—	—	57	3
Hemopexin	20,000	20,000	—	—	10 (stool)	100	100
Hepcidin	16	2	—	—	—	7	11
hCG alpha	4	4	—	—	—	12	12
HGF	2	2	—	—	—	12	12
HSP27/HspB1	4	4	—	—	—	12	12
HSP60	Neat	Neat	—	—	—	57	3
HSP70	—	2	—	—	—	12	12
HVEM/TNFRSF14	2	2	—	—	—	12	12
Hyaluordinase-1	32	32	—	—	—	57	3
IA-2/PTPRN	Neat	Neat	—	—	—	12	11
ICAM-1	1,000	1,000	—	—	—	101	101
ICAM-3	2	2	2	2	—	10	3
ICOS	Neat	Neat	—	—	—	57	3
ICOS-L/B7-H2	16	16	—	—	—	57	3
IgE	500	500	—	—	—	100	100
IGF-1	—	Neat	—	—	—	12	11
IGF-2R/CI-M6P-R	Neat	Neat	—	—	—	57	3
IGFBP-1	2	2	—	—	—	13	11
IGFBP-2	50	50	—	—	—	57	3
IGFBP-3	20	20	—	—	—	57	3
IGFBP-4	8	5	—	—	—	12	12
IGFBP-8	4	4	—	—	—	57	3
Ikaros	—	—	—	—	0.62 µg/well (cell lysate)	39 (with additives)	MSD Tris Wash Buffer (with additives)
IL-1R2	Neat	Neat	—	—	—	57	3
IL1RL1/ST2	4	4	—	—	—	7	37
IL-6R	50	50	50	50	—	7	8
IL-11	—	—	—	—	4 (Synovial Fluid)	101	3
IL-17RB	Neat	Neat	—	—	—	57	3
IL-18BP	4	4	—	—	—	57	3



Dash (—) indicates that the information for that sample type is not available.

18199-v15-2023Jun | 5

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES

## R-PLEX Antibody Set Dilution Factors and Assay/Antibody Diluent Combinations

Human							
Antibody Set	Dilution Factor					Primary Assay Diluent	Primary Antibody Diluent
	Serum	Plasma	Urine	CSF	Other		
IL-19	2	2	—	—	—	57	3
IL-22Rα2	Neat	Neat	—	—	—	57	3
IL-23R	Neat	Neat	—	—	—	57	3
IL-28B/IFN-A3	Neat	Neat	—	—	—	57	3
IL-32	Neat	Neat	—	—	—	12	11
IL-34	Neat	Neat	—	—	—	57	3
IL-36γ	Neat	Neat	—	—	—	13	11
IL-36RA	Neat	Neat	—	—	—	57	3
IL-37	Neat	Neat	—	—	—	57	3
Influenza A NP	—	—	—	—	2 (cell culture supernatant)	11	11
Influenza B NP	—	—	—	—	2 (cell culture supernatant)	11	11
IR (soluble)	4	4	—	—	—	100	100
IRS-1	—	—	—	—	0.63 µg/well (cell lysate)	13	12
Jagged-1	Neat	Neat	—	—	—	12	11
Kallikrein 3/PSA	Neat	Neat	—	—	—	57	3
Klotho	Neat	Neat	—	—	—	57	3
Lactotransferrin	1,000	1,000	—	—	—	100	100
LBP	500	500	—	—	—	12	11
LDL Receptor	4	4	—	—	—	12	11
LIF2	Neat	Neat	—	—	—	13	11
LIGHT/TNFSF14	Neat	Neat	—	—	—	57	3
LOX-1/OLR1	Neat	Neat	—	—	—	57	3
LOXL2	Neat	Neat	—	—	—	57	3
LRRK2	—	—	—	—	2.5–5.0 µg/well (cell lysate)	MSD Tris Lysis Buffer (with additives)	MSD Tris Wash Buffer (with additives)
LRRK2 (pS935)	—	—	—	—	2.5–5.0 µg/well (cell lysate)	MSD Tris Lysis Buffer (with additives)	MSD Tris Wash Buffer (with additives)
Lymphotactin	4	4	—	—	—	12	11
LYVE-1	100	100	—	—	—	57	3
MAdCAM-1	16	16	—	—	—	13	11
MCAM/CD146	Neat	Neat	—	—	—	57	3
Mcl-1/BAK Complex	—	—	—	—	5.0 µg/well (cell lysate)	39 (with additives)	39
Mcl-1/Bcl2-L-3	—	—	—	—	2.0 µg/well (cell lysate)	39 (with additives)	39
Mcl-1/BIM Complex	—	—	—	—	5.0 µg/well (cell lysate)	39 (with additives)	39
MerTK	Neat	Neat	—	—	—	57	3
Mesothelin	50	50	50	50	—	7	8
Met (soluble)	20	20	20	20	—	58	3
MIC-A	2	2	—	—	—	57	3
MIG	4	4	4	4	—	58	3
Midkine	2	2	—	—	—	13	11
MIP-4	1,000	1,000	—	—	—	57	3
MMP-1	10	10	10	10	—	7	3
MMP-2	2	2	—	—	—	58	3
MMP-3	10	10	10	10	—	7	3



Dash (—) indicates that the information for that sample type is not available.

18199-v15-2023Jun | 6

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES

## R-PLEX Antibody Set Dilution Factors and Assay/Antibody Diluent Combinations

Human							
Antibody Set	Dilution Factor					Primary Assay Diluent	Primary Antibody Diluent
	Serum	Plasma	Urine	CSF	Other		
MMP-7	2	2	—	—	—	12	12
MMP-8	2	2	—	—	—	58	3
MMP-9 (active)	16	8	—	—	—	57	3
MMP-9 (total)	40	40	—	—	—	58	3
MMP-10 (total)	Neat	Neat	—	—	—	12	11
MMP-12	Neat	Neat	—	—	—	2	3
MMP-13	Neat	Neat	—	—	—	57	3
MPIF-1/CCL23	2	2	—	—	—	57	3
MPO	20	20	—	—	—	57	3
MSP	50	50	—	—	—	57	3
MyD88	Neat	Neat	—	—	—	13	11
Myoglobin	50	50	—	—	—	13	11
Myosin BP (cardiac)	Neat	Neat	—	—	—	12	11
NAP-2/CXCL7	500	100	—	—	—	57	3
N-Cadherin	2	2	—	—	—	12	11
NCAM-1	100	100	—	—	—	101	3
Nectin-4	2	2	2	2	—	10	3
Nephrin	Neat	Neat	—	—	—	57	3
Neprilysin	2	2	—	—	—	12	11
Nesfatin-1	Neat	Neat	—	—	—	12	11
Neurofilament H	—	—	—	2	—	12	11
Neurofilament L	2	2	—	2	—	12	11
Neuronal pentraxin-1	10	10	—	100	—	12	11
Neuropilin-1	100	100	—	—	—	12	11
Neurotrophin-3	—	—	—	Neat	—	58	3
NGAL/LCN2	60	60	—	—	—	57	3
NPY	8	32	—	—	—	58	3
NT-ProBNP	50	50	—	—	—	58	3
Osteoactivin	10	10	10	10	—	7	3
Osteocalcin	50	50	50	50	—	7	8
Osteonectin	2	2	2	2	—	58	3
Osteopontin	—	4	—	—	—	7	37
Osteoprotegerin	2	2	2	2	—	58	3
OX40L	2	2	—	—	2 (cell lysate)	12	12
OXM/Glicentin	—	2	—	—	—	12	11
p38 (pT180/pY182)	—	—	—	—	20 µg/well (cell lysate)	39	39
p38 (total)	Neat	Neat	—	—	20 µg well (cell lysate)	39	39
PAF-AH	2	2	—	—	—	12	11
PC9/PCSK9	8	8	—	—	—	13	11
P-Cadherin	10	10	10	10	—	7	3
PDGF-A	2	16	—	—	—	57	3
PDGF-A/B	2	2	—	—	—	13	11
PDGF-B	2	2	—	—	—	7	11
PDGF-C	2	2	—	—	—	13	11
PDGFR-β (total)	8	5	—	—	—	101	101
Pentraxin 3	2	2	—	—	—	12	12



Dash (—) indicates that the information for that sample type is not available.

18199-v15-2023Jun | 7

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES

## R-PLEX Antibody Set Dilution Factors and Assay/Antibody Diluent Combinations

Human							
Antibody Set	Dilution Factor					Primary Assay Diluent	Primary Antibody Diluent
	Serum	Plasma	Urine	CSF	Other		
Perforin	8	8	—	—	—	58	3
Periostin/OSF-2	8	8	—	—	—	57	3
PF-4/CXCL4	500	100	—	—	—	57	3
PR/PGR	—	—	—	—	—	12	11
PRAS40 (total)	Neat	Neat	—	—	0.5 µg/well (cell lysate)	39	39
PRDX-1	—	4	—	—	—	7	37
PRDX-6	—	5	—	—	—	7	7
Presenilin-1 NTF	Neat	Neat	—	—	—	12	11
Procalcitonin	Neat	Neat	—	—	—	12	11
proCollagen alpha-1	8	8	—	—	—	12	11
Progranulin	4	4	—	—	—	57	3
Prolactin	10	10	—	—	—	13	12
proMMP-9	40	40	—	—	—	58	3
proMMP-10	2	—	—	—	—	13	11
proMMP-13	Neat	Neat	—	—	—	13	11
Protein C Receptor	Neat	Neat	—	—	—	12	11
P-Selectin	2	2	2	2	—	10	3
PSGL-1	—	4	—	—	—	7	37
PSP-D	4	4	—	—	—	57	3
PTEN (total)	Neat	Neat	—	—	20 µg/well (cell lysate)	39	39
PYY (active)	Neat	Neat	—	—	—	12	11
RAGE (soluble)	—	2	—	—	—	12	12
RANTES	50	50	50	50	—	7	8
RBP4	2,000	2,000	2	—	—	37	37
Reg-3-α	10	10	—	—	—	12	11
Relaxin-2	Neat	Neat	—	—	—	57	3
Renin	Neat	Neat	—	—	—	12	11
Resistin	50	50	50	50	—	58	3
S100A8/MRP8	2	2	—	—	—	12	12
S100A12	80	2	—	—	—	12	12
S100B	Neat	Neat	—	—	—	12	12
SAA	1,000	1,000	—	—	—	101	101
SARS-CoV-2 N	—	—	—	—	Neat (upper resp. sample)	11 (with additives)	11 (with additives)
SARS-CoV-2 Spike	—	—	—	—	Neat (upper resp. sample)	11 (with additives)	11 (with additives)
SCF	2	2	2	2	—	10	3
SCFR/Kit	50	50	—	—	—	7	11
Sclerostin	Neat	Neat	—	—	—	13	11
Serpin A4	100,000	100,000	—	—	—	57	3
Serpin A12/Vaspin	2	2	—	—	—	12	12
Serpin E1 (active)	16	16	—	—	—	57	3
Serpin E1 (inactive)	30	30	—	—	—	57	3
Serpin E1 (total)	30	30	—	—	—	57	3
Serpin F1/PEDF	4	4	—	—	—	101	3
Serpin F2	20,000	20,000	—	—	2 (stool)	100	100
SHBG	1,000	—	—	—	—	57	3



Dash (—) indicates that the information for that sample type is not available.



## R-PLEX Antibody Set Dilution Factors and Assay/Antibody Diluent Combinations

Human							
Antibody Set	Dilution Factor					Primary Assay Diluent	Primary Antibody Diluent
	Serum	Plasma	Urine	CSF	Other		
SOD2	Neat	Neat	—	—	20 µg/well (cell lysate)	39	39
SPINK-1	2	2	—	—	—	57	3
SPK 1	Neat	Neat	—	—	—	57	3
SR-PSOX/CXCL16	50	50	—	—	—	12	11
sTfR-1	—	5,000	—	—	—	101	37
Syndecan-1	2	2	—	—	—	57	3
Syndecan-4	4	4	—	—	—	57	3
TACI/TNFRSF13	8	8	—	—	—	57	3
Tau (total)	—	—	—	4	—	35	35
TECK/CCL25	Neat	Neat	—	—	—	58	3
Tenascin C	4,000	4,000	4,000	4,000	—	12	12
TFF2	8	8	—	—	—	12	11
TFF3	—	4	—	—	—	7	37
TFPI	4	4	—	—	—	12	11
TGF-α	Neat	Neat	—	—	—	12	11
TGF-β RII	Neat	Neat	—	—	—	12	11
TGM-2	2	2	—	—	—	12	11
Thrombomodulin	2	2	2	2	—	10	3
Thyroglobulin	2	2	—	—	—	12	12
TIMP-1	—	100	—	—	—	101	37
TIMP-3	Neat	Neat	—	—	—	12	11
TIMP-4	4	4	—	—	—	12	11
TNFAIP6/TSG-6	2	2	—	—	—	12	12
TNF-RI	10	10	10	10	—	7	3
TNF-RII	10	10	—	—	—	58	3
TNFRSF10C	20	20	—	—	—	12	12
tPA	4	4	—	—	—	57	3
Troponin I (cardiac)	4	4	—	—	—	7	37
Troponin T (cardiac)	20	—	—	—	—	7	22
Trk-A (total)	Neat	Neat	—	—	20 µg/well (cell lysate)	39	39
TSP-1	2	2	—	—	—	12	37
TSP-2	8	8	—	—	—	12	11
TWEAK/TNFSF12	Neat	Neat	—	—	—	13	11
TYRO-3	4	4	—	—	—	12	11
U-PAR	—	2	—	—	—	12	12
Uromodulin	500	500	500	—	—	37	37
VCAM-1	1,000	1,000	—	—	—	101	101
VE-Cadherin	100	100	—	—	—	12	11
VEGFR-2/KDR	50	50	—	—	—	7	11
VEGFR-3/Flt-4	2	2	—	—	—	57	3
VILIP-1	—	2	—	—	—	101	101
Visfatin	2	2	—	—	—	35	35
VISTA/B7-H5	Neat	Neat	—	—	—	57	3
vWF	—	1,000	—	—	—	12	12
ZnT8	Neat	Neat	—	—	—	12	11



Dash (—) indicates that the information for that sample type is not available.

18199-v15-2023Jun | 9

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES

## R-PLEX Antibody Set Dilution Factors and Assay/Antibody Diluent Combinations

Nonhuman Primate							
Antibody Set	Dilution Factor					Primary Sample Diluent	Primary Antibody Diluent
	Serum	Plasma	Urine	CSF	Other		
Alpha-amylase 2A	5,000	5,000	—	—	—	58	3
CD177	16	—	—	—	—	58	3



Dash (—) indicates that the information for that sample type is not available.

## R-PLEX Antibody Set Dilution Factors and Assay/Antibody Diluent Combinations

Mouse							
Antibody Set	Dilution Factor					Primary Sample Diluent	Primary Antibody Diluent
	Serum	Plasma	Urine	CSF	Other		
Adiponectin	1,000	1,000	—	—	—	100	100
Amphiregulin	Neat	Neat	—	—	—	41	45
AXL/UFO	16	16	—	—	—	41	45
C1q receptor	500	500	—	—	—	41	45
CCL6	Neat	16	—	—	—	41	45
CD14	2	2	—	—	—	41	45
CD163	Neat	Neat	—	—	—	13	3
Clusterin	1,000	1,000	—	—	—	41	45
Complement C5a	10,000	8	—	—	—	41	45
CRP	1000	1000	—	—	—	41	45
CTLA-4	Neat	Neat	—	—	—	13	3
Cystatin C	500	500	—	—	—	100	100
Elastase-2	2	Neat	—	—	—	41	45
Eotaxin-2	Neat	Neat	—	—	—	41	45
Fractalkine	2	2	—	—	—	41	45
FTL3L	Neat	Neat	—	—	—	13	3
GAS-6/AXL-L	64	64	—	—	—	41	45
Galectin-3	64	64	—	—	—	42	40
G-CSF	2	2	—	—	—	57	3
GDF-15	Neat	Neat	—	—	—	13	3
Granzyme B	Neat	Neat	—	—	—	13	11
HAVCR1/KIM-1	Neat	Neat	—	—	—	41	45
HGF	2	2	—	—	—	41	45
ICAM-1	50	50	—	—	—	41	45
IGF-1	256	64	—	—	—	41	45
IGFBP-5	16	16	—	—	—	13	3
IL-1 $\alpha$	Neat	Neat	—	—	—	57	3
IL-1RA	Neat	Neat	—	—	—	12	11
IL-6R $\alpha$	16	16	—	—	—	41	45
IL-7	Neat	Neat	—	—	—	57	3
IL-18	Neat	Neat	—	—	—	42	40
IL-24	Neat	Neat	—	—	—	13	3
I-TAC	Neat	Neat	—	—	—	13	3
LIX/CXCL5	Neat	Neat	—	—	—	41	45
M-CSF	16	16	—	—	—	57	3
MIF	Neat	Neat	—	—	—	42	40
MIG	2	2	—	—	—	41	45
MIP-1 $\gamma$ /CCL9	256	256	—	—	—	41	45
MMP-3 (total)	2	2	—	—	—	41	45
Oncostatin-M	Neat	Neat	—	—	—	41	45
Osteoprotegerin	4	4	—	—	—	41	45
PDGF-B	4	Neat	—	—	—	13	3
PD-L1	Neat	Neat	—	—	—	13	3
Pentraxin 3	8	8	—	—	—	42	40
Periostin/OSF-2	100	100	—	—	—	41	45
PF-4/Cxcl4	5,000	2,500	—	—	—	41	45
Prolactin	Neat	Neat	—	—	—	13	3



Dash (—) indicates that the information for that sample type is not available.

18199-v15-2023Jun | 11

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES

## R-PLEX Antibody Set Dilution Factors and Assay/Antibody Diluent Combinations

Mouse							
Antibody Set	Dilution Factor					Primary Sample Diluent	Primary Antibody Diluent
	Serum	Plasma	Urine	CSF	Other		
<b>PSP-D</b>	2	2	—	—	—	41	45
<b>RBP4</b>	1,000	1,000	—	—	—	41	45
<b>Resistin</b>	20	20	—	—	—	11	11
<b>S100A8/MRP8</b>	—	Neat	—	—	—	42	40
<b>SAA</b>	500	500	—	—	—	41	45
<b>Tau (total)</b>	—	—	—	4	—	35	35
<b>TIMP-1</b>	Neat	Neat	—	—	—	57	3
<b>TNF-RII</b>	32	32	—	—	—	100	100
<b>TREM-1</b>	2	2	—	—	—	41	45
<b>VCAM-1</b>	1,000	1,000	—	—	—	41	45
<b>VEGFR-1/Flt-1</b>	Neat	Neat	—	—	—	12	11
<b>YM1</b>	16	4	—	—	—	41	45



Dash (—) indicates that the information for that sample type is not available.

## R-PLEX Antibody Set Dilution Factors and Assay/Antibody Diluent Combinations

Rat							
Antibody Set	Dilution Factor					Primary Sample Diluent	Primary Antibody Diluent
	Serum	Plasma	Urine	CSF	Other		
A2M	20,000	20,000	—	—	—	100	100
Adiponectin	1,000	1,000	—	—	—	100	100
AGP	20,000	20,000	—	—	—	100	100
Albumin	—	—	8	—	—	29	29
β-NGF	Neat	Neat	—	—	—	41	45
Cystatin C	500	500	—	—	—	100	100
EPO	Neat	Neat	—	—	—	42	40
Fractalkine	Neat	Neat	—	—	—	42	40
GM-CSF	2	2	—	—	—	42	40
HAVCR1/KIM-1	4	4	—	—	—	42	40
IFN-γ	4	4	4	—	—	42	40
IL-1α	4	4	—	—	—	42	40
IL-1β	4	4	—	—	—	42	40
IL-2	2	2	—	—	—	42	40
IL-4	4	4	—	—	—	42	40
IL-5	2	2	—	—	—	42	40
IL-6	4	4	—	—	—	42	40
IL-10	2	2	—	—	—	42	40
IL-13	Neat	Neat	—	—	—	42	40
KC/GRO	8	8	—	—	—	42	40
MCP-1	Neat	Neat	—	—	—	42	40
MIP-3α	Neat	Neat	—	—	—	42	40
NGAL/LCN2	100	100	—	—	—	7	100
Osteopontin	—	—	10	—	—	29	100
Resistin	20	20	—	—	—	11	11
TIMP-1	20	20	—	—	—	11	5
TNF-α	2	2	—	—	—	42	40
TNF-RII	32	32	—	—	—	100	100
TREM-1	2	2	—	—	—	41	45
VEGF-A	Neat	Neat	—	—	—	42	40

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. ©2015-2018, 2020-2023 Meso Scale Diagnostics, LLC. All rights reserved.



Dash (—) indicates that the information for that sample type is not available.

18199-v15-2023Jun | 13  
FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES