

Quanterix SR-X™ Ultrasensitive Biomarker Assay Menu

Singleplex Advantage Assays

Analyte	LoD (pg/mL)	LoQ (pg/mL)	Dynamic Range (pg/mL)	Median Endogenous (pg/mL)	Sample Volume	Sample Type [†]	Catalog Number
Aβ40	0.170	1.24	0-800	177	25 µL	C, E	101672
Aβ42	0.035	0.137	0-400	10.2	25 µL	C, E	101664
C-Peptide	0.008	0.027	0-400	1.53 (Type I Diabetes)	25 µL	C, E	100199
HIV p24	0.004	0.016	0-30	N/A	125 µL	E, S	102215
IFN-γ	0.014	0.049	0-800	0.33	65 µL	E, S	103337
IL-5	0.0034	0.0165	0 - 12	0.210	76 µL	E, S	102860
IL-6	0.006	0.010	0-120	1.27	25 µL	E, S	101622
IL-8	0.038	0.146	0-300	4.27	25 µL	E, S	100198
IL-10	0.005	0.041	0-120	1.29	25 µL	E, S	101643
IL-13	0.001	0.007	0-30	0.069	50 µL	E, S	102732
IL-15	0.002	0.007	0-40	3.23	25 µL	E, S	100794
IL-17A	0.011	0.011	0-120	0.073	25 µL	E, S	101599
NF-light™ for SR-X	0.072	0.343	0-1,800	7.34	25 µL	C, E, S	103400
PSA	0.006	0.028	0-400	1.81	25 µL	E, S	101478
pTau-231	0.284	1.23	0-1,200	47.5	25 µL	C	102292
pTau-181	0.756	1.20	0-85	33.1	25 µL	C	103377
pTau-181 V2	0.041	0.085	0-412	0.990	25 µL	C, E, S	103714
SNAP-25	1.06	3.20	1-1,000	59.6	32.5 µL	C	103575
Tau	0.014	0.049	0-400	2.65	38 µL	C, E, S	101552
TDP43	0.780	8.23	0-8,000	191	25 µL	C, E, S	103293
TNF-α	0.003	0.017	0-200	1.48	25 µL	E, S	101580

Singleplex Discovery Assays

Analyte	LoD (pg/mL)	LoQ (pg/mL)	Dynamic Range (pg/mL)	Median Endogenous (pg/mL)	Sample Volume	Sample Type [†]	Catalog Number
α-Synuclein	0.440	4.12	0-10,000	3,976	15.2 µL	C, E, S	102233
BDNF	0.021	0.068	0-6,400	4,578	5 µL	C, E, S	102039
GFAP	0.26	1.37	0-4,000	68.4	38 µL	C, E, S	102336
IL-22 (total)	0.004	0.021	0-120	6.62	25 µL	E, S	103071
MMP-9	1.09	4.88	0-5,000 ng/ml	680	5 µL	C, P, S	102491
Mouse IL-1β	0.0081	0.0309	0-120	0.296	25 µL	E, S	102517
Mouse IL-6	0.072	0.240	0-700	38.34	25 µL	E, S	102869
Mouse Tau	0.428	0.82	0-2,400	19.6	25 µL	C, E, S	102209
PD-L1	0.048	0.052	0-4,300	40.1	5 µL	E, S	102648

[†]C = CSF, E = EDTA plasma, P=platelet poor heparin plasma, S = serum

Multiplex Advantage Assays - Human

Assay	Analytes	LoD (pg/mL)	LoQ (pg/mL)	Dynamic Range (pg/mL)	Median Endogenous (pg/mL)	Sample Volume	Sample Type [†]	Catalog Number
Neurology 2-Plex B	GFAP	0.410	4.15	0-40,000	43.2	25 µL	C, E, S	103520
	NF-light	0.071	0.40	0-2000	3.70	5 µL (CSF)		
Neurology 3-Plex A	Aβ40	0.196	0.675	0-800	209	25 µL 5 µL (CSF)	C, E	101995
	Aβ42	0.045	0.142	0-400	11.1			
	Tau	0.019	0.063	0-400	1.43			
Neurology 4-Plex A	GFAP	0.221	0.467	0-4,000	89.7	38 µL 5 µL (CSF)	C, E, S	102153
	NF-light	0.104	0.241	0-2000	10.6			
	Tau	0.024	0.053	0-400	2.21			
	UCH-L1	1.74	5.45	0-40,000	12.21			
Neurology 4-Plex B	GFAP	1.51	9.38	0-40,000	50.0	25 µL 5 µL (CSF)	C, E, S	103345
	NF-light	0.105	0.500	0-2,000	5.91			
	Tau	0.041	0.125	0-400	0.95			
	UCH-L1	1.90	9.38	0-40,000	6.81			

[†]C = CSF, E = EDTA plasma, S = serum

Simoa Advantage assay performance and precision specifications are determined according to Clinical Laboratory Standards Institute guidelines for consistency, sensitivity, precision, and robustness. Details are included in each kit's product insert and certificate of analysis. Kit includes calibrator concentrates, calibrator diluents, beads, detectors, SBG, RGP, and sample diluents. Validation testing includes:

- Dynamic range
- Sensitivity
- Calibration curve fitting
- Intra-run, inter-plate, and inter-lot precision and accuracy
- Robustness and stability of calibrators, antibodies, and controls
- Analysis of normal samples for the specified species to determine the normal range of biomarker concentration detected with the assay
- Specificity, spike recovery, and dilution linearity

Simoa Discovery assay kits are provided as complete kits packaged with reagents for processing up to 192 samples. Includes calibrator concentrate, calibrator diluent, bead stock, bead diluent, detector stock, detector diluent, SBG concentrate, SBG diluent, RGP, and sample diluent. Representative performance is determined in biological matrices and includes:

- Dynamic range
- Sensitivity
- Calibration curve fitting
- Intra-run, inter-plate, and inter-lot precision and accuracy
- Robustness and stability of calibrators, antibodies, and controls
- Analysis of normal samples for the specified species to determine the normal range of biomarker concentration detected with the assay
- Specificity, spike recovery, and dilution linearity