

Product Number:	102153
Lot Number:	501960
Expiration:	23-Feb-2020
Platform(s):	HD-1/HD-X & SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102141	919210
Detector Reagent	102143	919215
SBG Reagent	102145	905402
Sample Diluent	102147	917503
Calibrators	102150	920401
Controls	102151-102152	920402
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ Neuro 4-Plex A SR-X Kit Instructions (KI-0031).

Release Materials

Tau

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920402	1.58	1.14-1.71
Control 2	920402	65.7	56.6-85.0
Panel 1	MF-101617-01	1.11	0.605-1.92
Panel 2	MF-101617-02	25.8	22.7-44.5
Panel 3	MF-101617-03	121	96.8-194

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920402	4.70	4.25-6.38
Control 2	920402	426	368-552
Panel 1	MF-101617-01	4.55	3.08-6.38
Panel 2	MF-101617-02	36.3	16.7-50.4
Panel 3	MF-101617-03	366	234-541

GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920402	19.3	16.6-24.9
Control 2	920402	930	799-1199
Panel 1	MF-101617-01	53.1	40.9-62.1
Panel 2	MF-101617-02	168	111-227
Panel 3	MF-101617-03	2166	2085-3303

UCHL1®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920402	111	78.9-140
Control 2	920402	9482	7735-11602
Panel 1	MF-101617-01	54.5	42.9-87.8
Panel 2	MF-101617-02	276	249-503
Panel 3	MF-101617-03	4548	4232-6348

Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Review/Approval Patrizia Stadler Assoc. Director, QC p.stadler 31 JUL 2019
Name Title Signature/Date