

Product Number:	102153
Lot Number:	502001
Expiration:	19-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	922801
Detector Reagent	102143	922803
SBG Reagent	102145	908107
Sample Diluent	102147	905009
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	1.66	1.10-1.71
Control 2	920402	70.6	57.1-85.6
Panel 1	MF-101617-01	1.36	0.605-1.92
Panel 2	MF-101617-02	30.5	22.7-44.5
Panel 3	MF-101617-03	128	96.8-194

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920402	5.43	4.23-6.34
Control 2	920402	402	371-556
Panel 1	MF-101617-01	5.10	3.08-6.38
Panel 2	MF-101617-02	36.5	16.7-50.4
Panel 3	MF-101617-03	364	234-541

GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920402	19.7	16.5-24.7
Control 2	920402	822	809-1213
Panel 1	MF-101617-01	45.4	40.9-62.1
Panel 2	MF-101617-02	185	111-227
Panel 3	MF-101617-03	2165	2085-3303

UCHL1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920402	89.3	78.3-138
Control 2	920402	8632	7808-11713
Panel 1	MF-101617-01	59.4	42.9-87.8
Panel 2	MF-101617-02	408	249-503
Panel 3	MF-101617-03	4673	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	<i>P. Stodler</i>	28 Aug 2019
	Name	Title	Signature/Date	