

Product Number:	103520
Lot Number:	502760
Expiration:	15-Feb-2022
Platform(s):	HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	103510	104612
Detector Reagent	103512	104609
SBG Reagent	103514	104707
Sample Diluent	103516	104711
Calibrators	103517	104902 (A-E), 106910 (F-H)
Control 1	103518	104903
Control 2	103519	104903
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 2-Plex B SR-X Kit Instructions (KI-0062).

Release Materials

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	104903	6.71	5.23-7.85
Control 2	104903	653	524-786
Panel 1	929401	3.83	2.63-4.73
Panel 2	929402	6.65	4.56-10.5
Panel 3	929403	86.2	74.3-112
Panel 4	929404	365	348-531

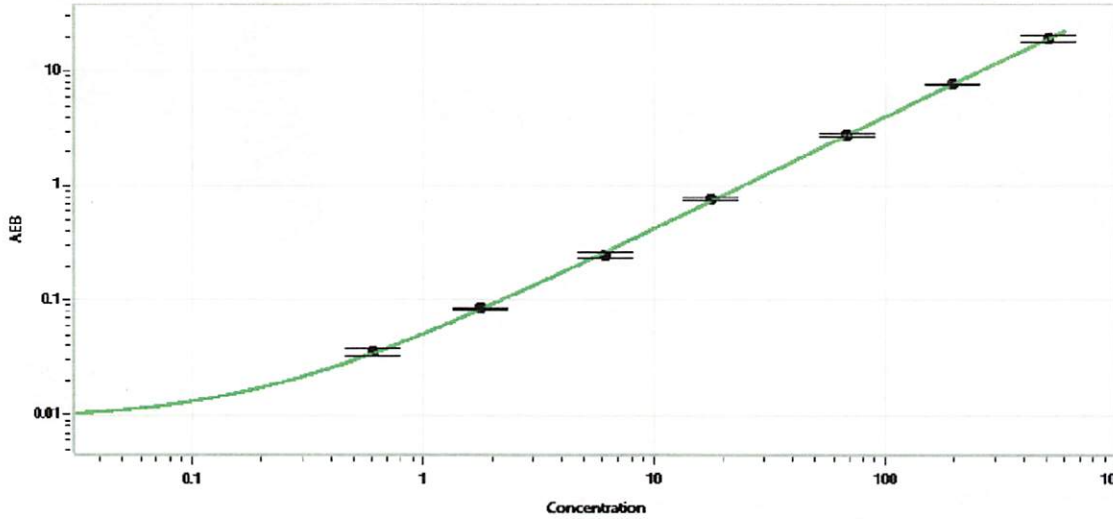
GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	104903	76.7	60.9-91.3
Control 2	104903	8873	6946-10419
Panel 1	929401	51.6	39.8-64.9
Panel 2	929402	45.8	34.1-71.7
Panel 3	929403	1189	1007-1511
Panel 4	929404	4048	3421-5189

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

Review/Approval	Junpil Park	QC Supervisor		07 May 2021
	Name	Title	Signature/Date	

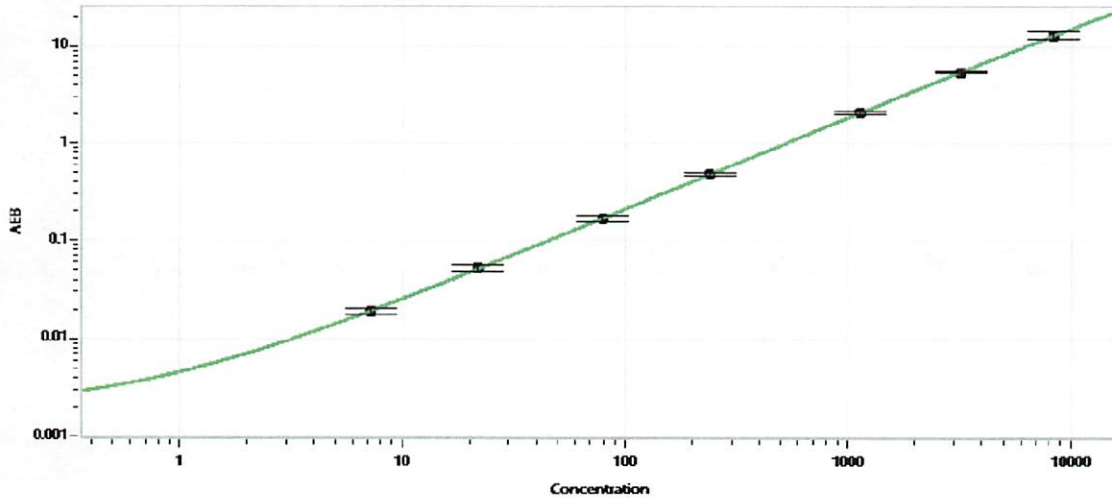
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.611
C	1.77
D	6.14
E	17.6
F	68.0
G	197
H	510

GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	7.19
C	21.8
D	79.6
E	239
F	1126
G	3230
H	8314