

Product Number:	103520
Lot Number:	503094
Expiration:	19-Jul-2022
Platform(s):	HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103510	123003
Detector Reagent	103512	124512
SBG Reagent	103514	123010
Sample Diluent	103516	123011
Calibrators	103517	113410
Control 1	103518	113411
Control 2	103519	113411
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).



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

Release Materials

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	113411	7.89	5.80-8.70
Control 2	113411	694	591-887
Panel 1	929401	3.59	2.63-4.73
Panel 2	929402	6.80	4.56-10.5
Panel 3	929403	84.1	74.3-112
Panel 4	929404	351	325-554

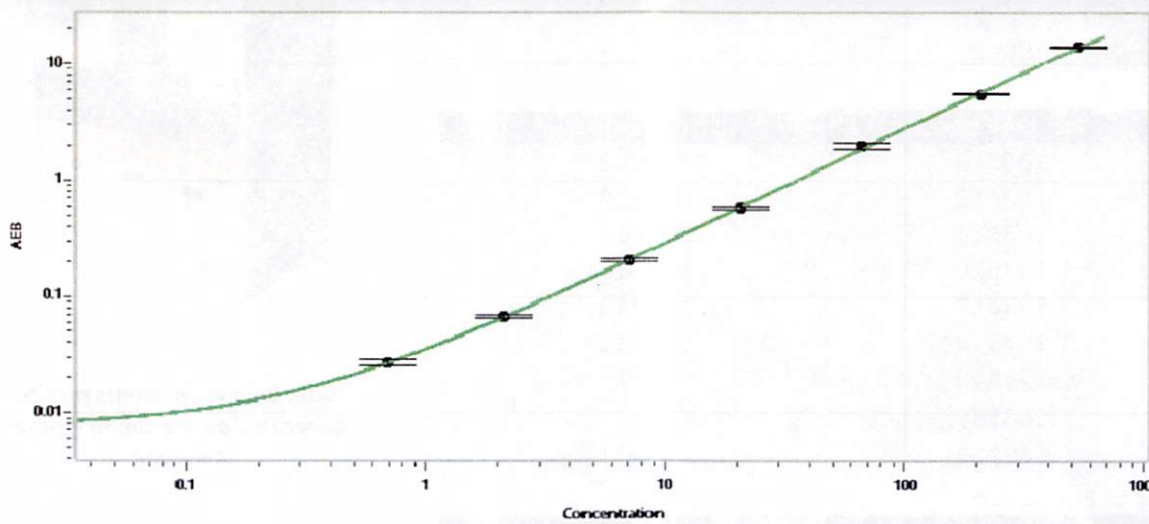
GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	113411	109	86.6-130
Control 2	113411	10623	9326-13990
Panel 1	929401	55.1	39.8-64.9
Panel 2	929402	53.8	34.1-71.7
Panel 3	929403	1236	1007-1511
Panel 4	929404	3669	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 [Signature] 17 Sep 2024
Name Title Signature/Date

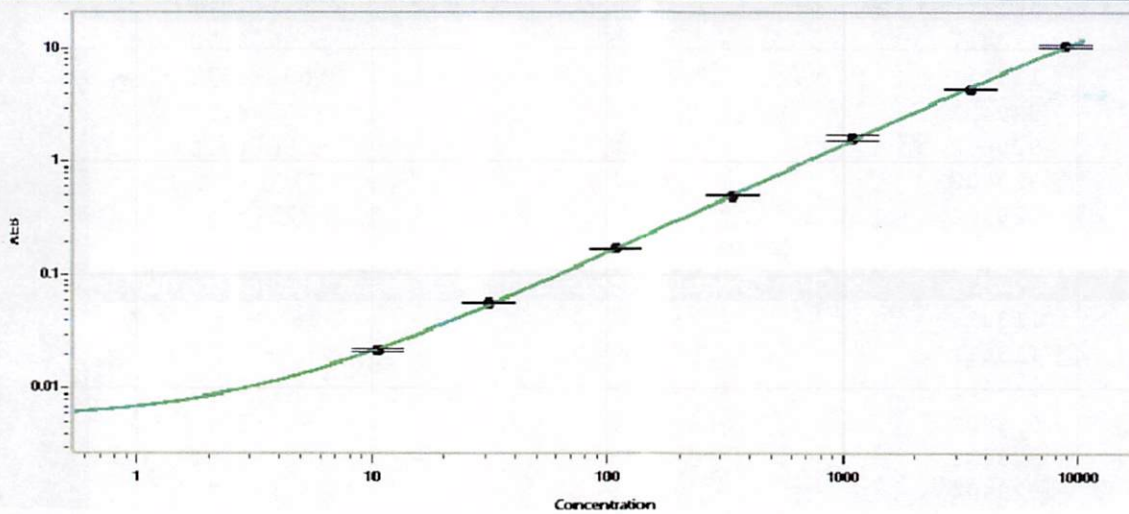
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.688
C	2.13
D	7.04
E	20.6
F	66.1
G	209
H	528

GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	10.6
C	31.2
D	108
E	339
F	1095
G	3480
H	8791