

Product Number:	103520
Lot Number:	502713
Expiration:	23-Nov-2021
Platform(s):	HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103510	102914
Detector Reagent	103512	103005
SBG Reagent	103514	103006
Sample Diluent	103516	034417
Calibrators	103517	008701
Control 1	103518	008702
Control 2	103519	008702
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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).



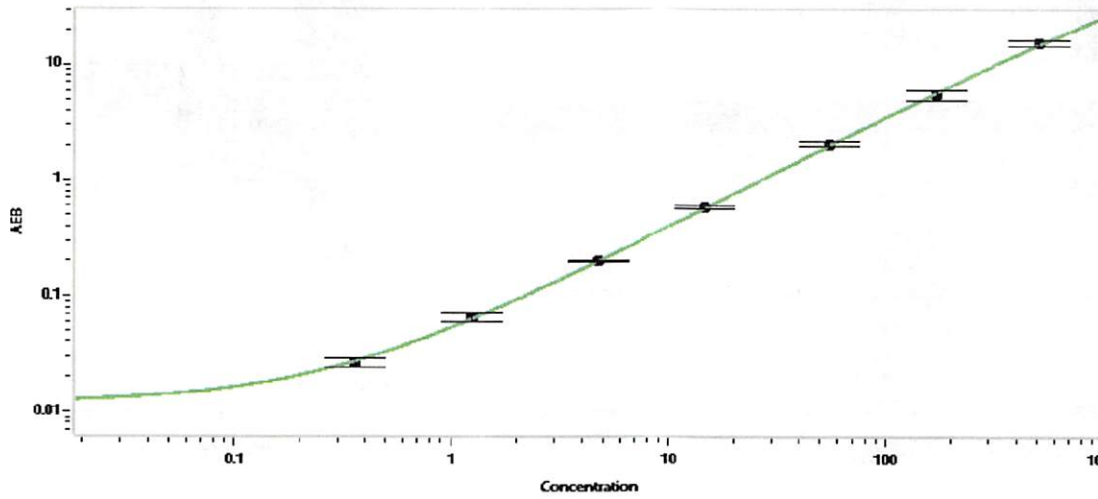
Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials			
NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	008702	5.39	3.45-7.62
Control 2	008702	611	482-851
Panel 1	929401	3.34	2.50-4.78
Panel 2	929402	6.75	4.72-10.8
Panel 3	929403	87.9	75.7-114
Panel 4	929404	428	367-550
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	008702	90.1	67.6-116
Control 2	008702	10035	8194-14792
Panel 1	929401	51.6	39.9-65.2
Panel 2	929402	49.1	36.9-72.7
Panel 3	929403	1251	1021-1532
Panel 4	929404	4255	3495-5243

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

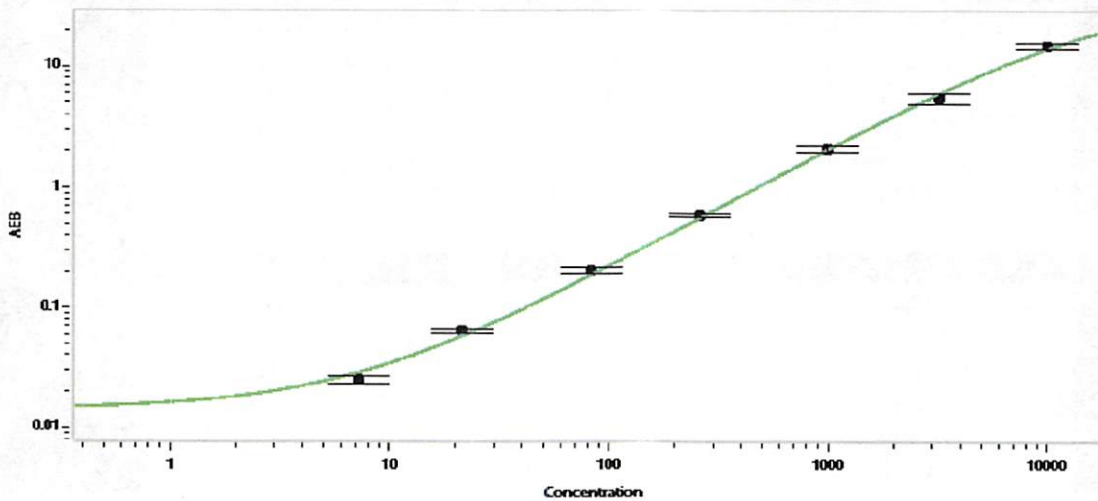
Review/Approval	Junpil Park	QC Supervisor		12 Feb 2021
	Name	Title	Signature/Date	

**NF-light® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	0.364
C	1.25
D	4.75
E	14.5
F	54.6
G	171
H	509

**GFAP® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	7.17
C	21.6
D	83.7
E	260
F	990
G	3175
H	9950