

Product Number:	103520
Lot Number:	503155
Expiration:	5-Aug-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	123003
Detector Reagent	103512	124512
SBG Reagent	103514	123010
Sample Diluent	103516	123011
Calibrators	103517	130401
Control 1	103518	130402
Control 2	103519	130402
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).

**Release Materials**

**NF-light®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	130402	8.81	6.97 - 10.5
Control 2	130402	828	669 - 1003
Panel 1	128902	13.1	9.26 - 13.9
Panel 2	128902	12.7	9.33 - 14.9
Panel 3	128902	138	101 - 151
Panel 4	128902	367	279 - 419

**GFAP®**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	130402	110	83.1 - 125
Control 2	130402	11815	9141 - 13712
Panel 1	128902	81.8	45.1 - 93.6
Panel 2	128902	70.6	38.3 - 91.2
Panel 3	128902	5186	3411 - 5538
Panel 4	128902	21623	16718 - 25077

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

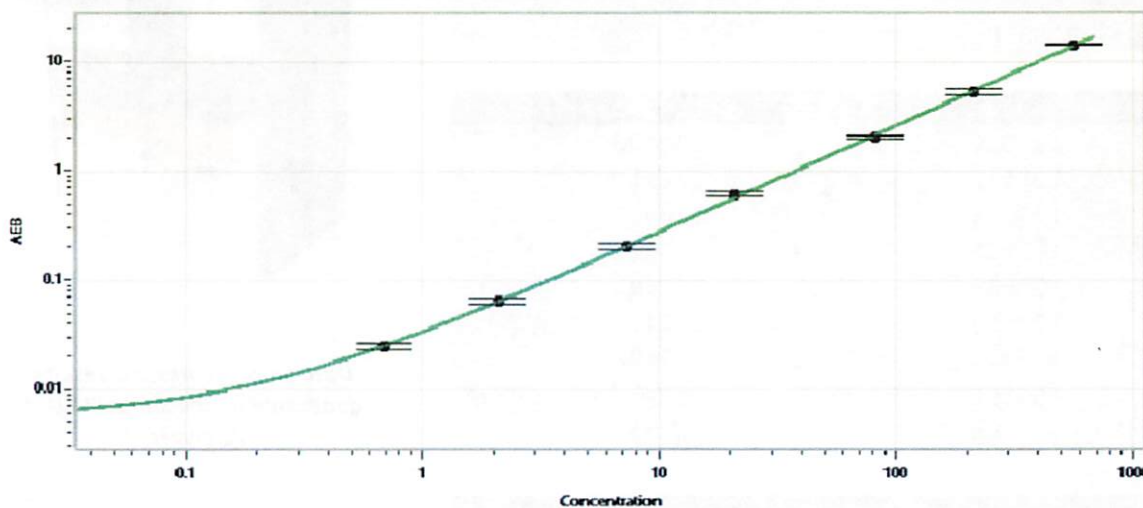
Review/Approval

Junpil Park  
Name

QC Supervisor  
Title

 10 NOV 2021  
Signature/Date

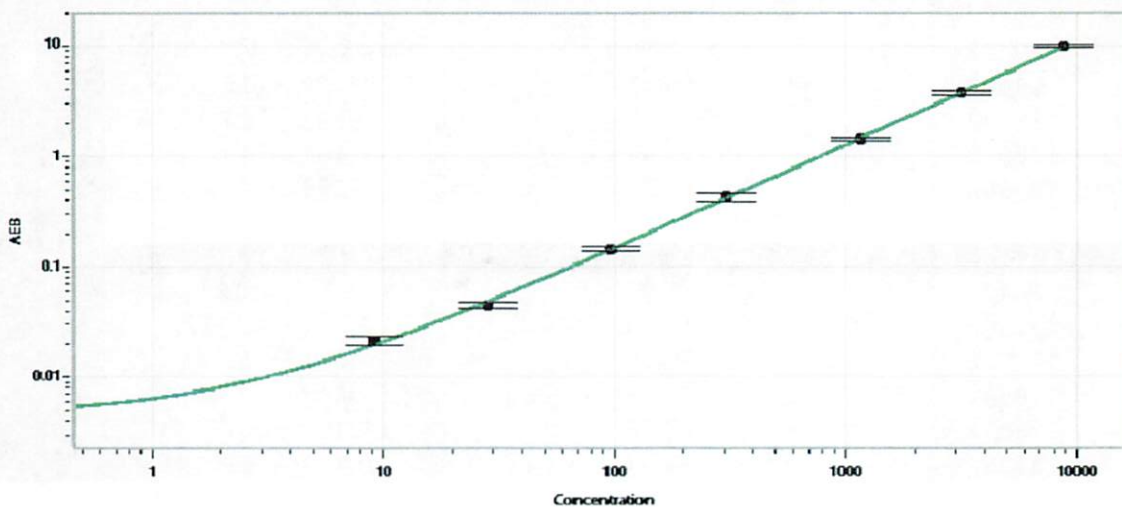
**NF-light® Calibration Curve**



Calibrator Levels  
(pg/mL)

A	0.000
B	0.687
C	2.09
D	7.31
E	20.9
F	82.0
G	214
H	563

**GFAP® Calibration Curve**



Calibrator Levels  
(pg/mL)

A	0.000
B	9.19
C	28.6
D	97.9
E	306
F	1183
G	3206
H	8794