

Product Number:	103520
Lot Number:	503233
Expiration:	18-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	134901
Control 1	103518	134902
Control 2	103519	134902
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	134902	4.37	3.32 - 5.39
Control 2	134902	405	335 - 503
Panel 1	128902	14.2	7.33 - 16.3
Panel 2	128902	12.3	8.91 - 15.3
Panel 3	128902	141	101 - 152
Panel 4	128902	362	261 - 436

GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134902	99.1	77.3 - 146
Control 2	134902	12491	9453 - 15447
Panel 1	128902	74.7	44.4 - 101
Panel 2	128902	68.5	45.0 - 79.6
Panel 3	128902	4924	3319 - 6109
Panel 4	128902	22328	17135 - 25703

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

Review/Approval

Wenkin Tsa.

Name

QC

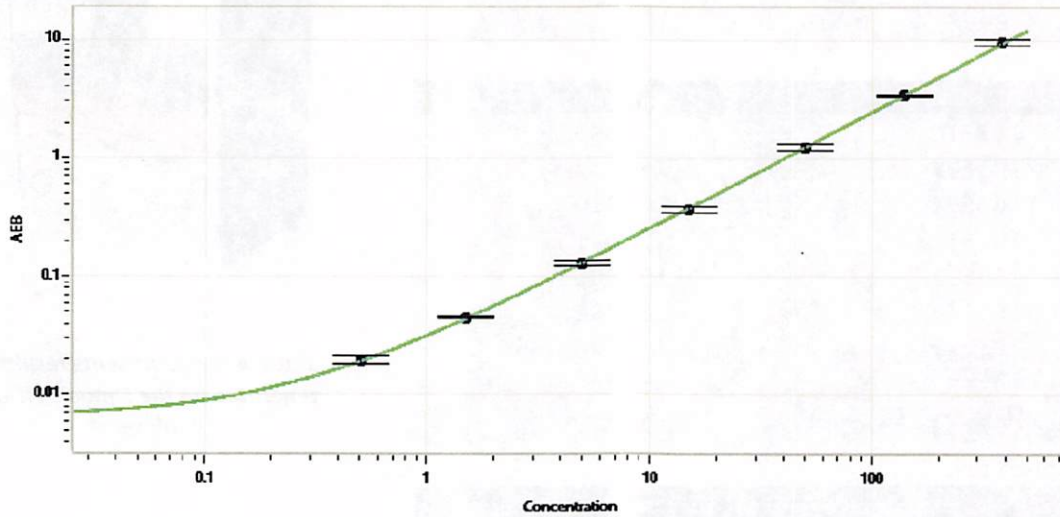
Title

[Handwritten Signature]

11 Jun 2022

Signature/Date

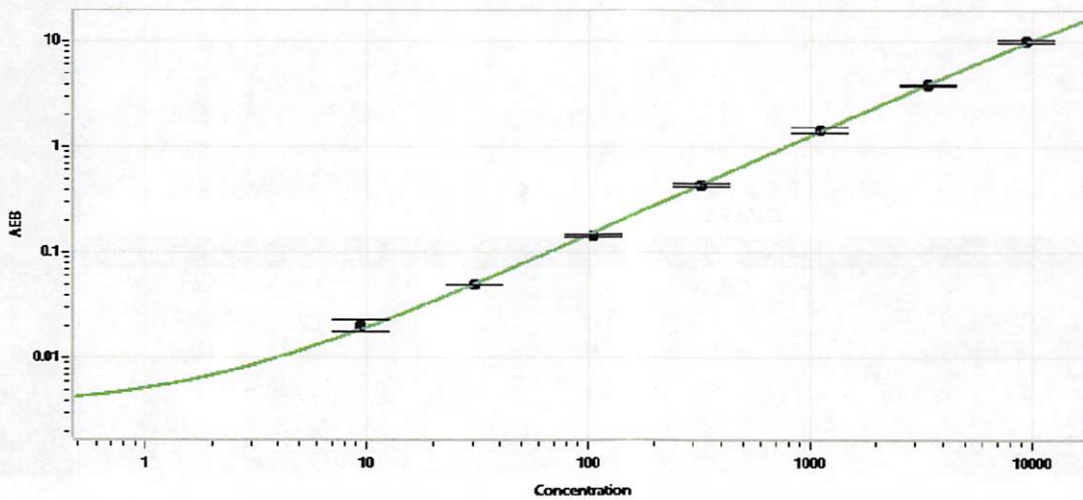
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.506
C	1.49
D	4.95
E	15.1
F	50.3
G	139
H	381

GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	9.40
C	30.9
D	105
E	323
F	1103
G	3397
H	9403