

Product Number:	103520
Lot Number:	502297
Expiration:	21-Jan-2021
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	006203
Detector Reagent	103512	006515
SBG Reagent	103514	006201
Sample Diluent	103516	002116
Calibrators	103517	008701
Controls	103518-103519	008702
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	008702	4.49	3.23-7.95
Control 2	008702	622	495-875
Panel 1	929401	3.17	2.56-4.94
Panel 2	929402	5.96	5.09-11.1
Panel 3	929403	90.5	77.5-116
Panel 4	929404	429	375-562

GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	008702	85.1	73.9-111
Control 2	008702	11077	9108-14850
Panel 1	929401	50.2	38.9-66.8
Panel 2	929402	44.1	39.3-74.2
Panel 3	929403	1244	1028-1542
Panel 4	929404	3976	3500-5314

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

Review/Approval

Shivani Goel

QC Manager

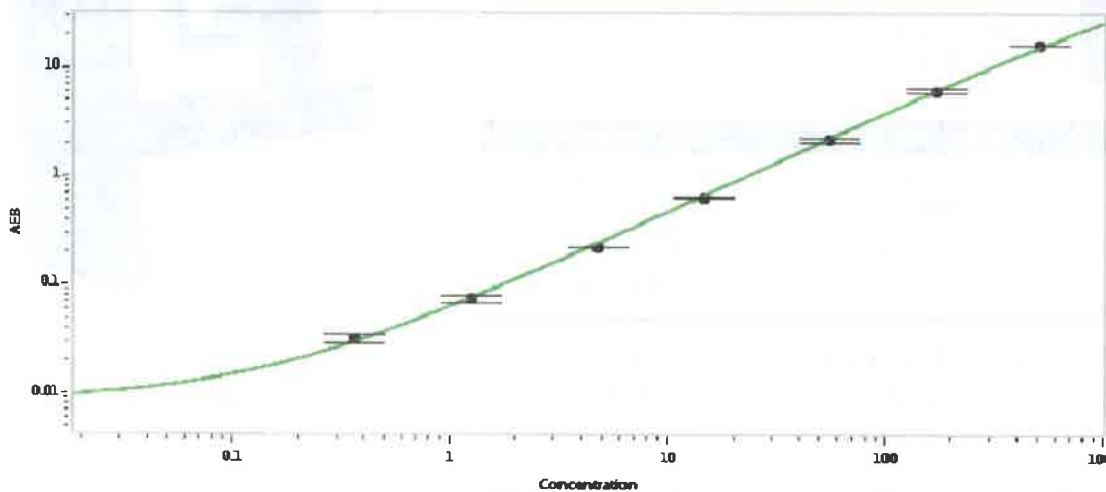
Shivani Goel 01 May 2020

Name

Title

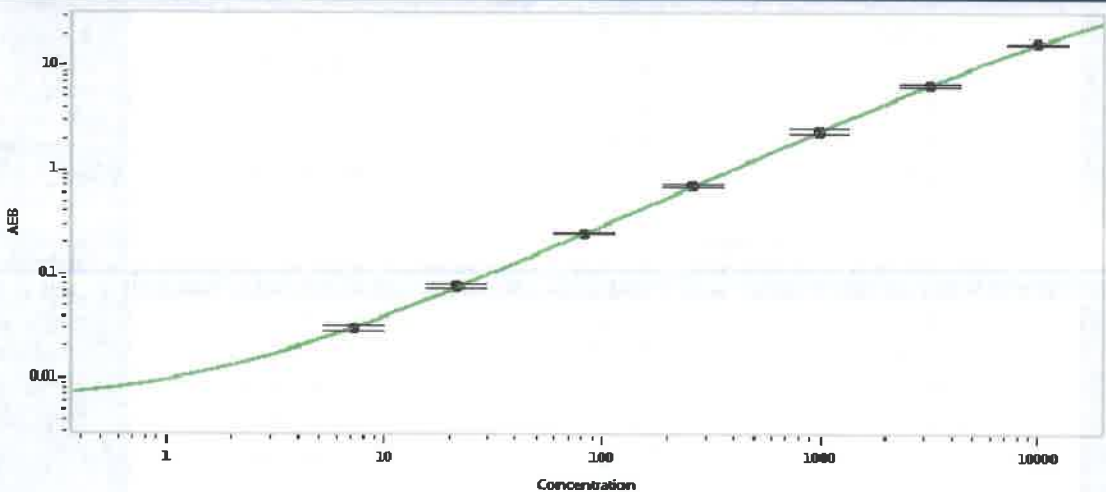
Signature/Date

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