

Product Number:	103520
Lot Number:	503866
Expiration:	4-Aug-2024
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	323611
Detector Reagent	103512	323702
SBG Reagent	103514	323703
Sample Diluent	103516	323612
Calibrators	103517	320202
Control 1	103518	320203
Control 2	103519	320203
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP / SBG Diluent (SR-X) are 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)	Spec. Range (pg/mL)
Control 1	320203	7.28	5.56 - 8.34
Control 2	320203	578	439 - 692
Panel 1	310007	7.67	4.60 - 9.38
Panel 2	323304	6.34	4.61 - 6.91
Panel 3	306904	88.3	74.0 - 111
Panel 4	306904	419	359 - 597

**GFAP®**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	320203	112	82.4 - 137
Control 2	320203	11394	8812 - 13219
Panel 1	310007	78.9	45.6 - 86.0
Panel 2	323304	62.1	31.9 - 72.1
Panel 3	306904	1078	795 - 1440
Panel 4	306904	7803	6199 - 9299

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

Review/Approval

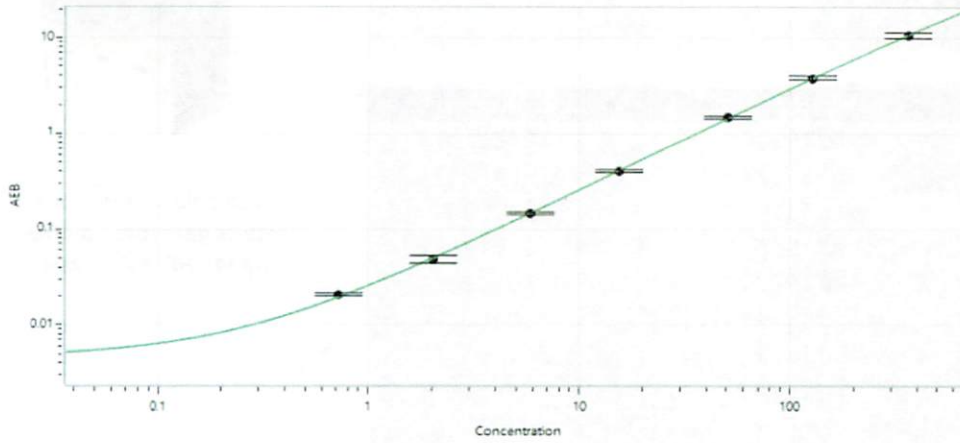
*Sonia Khillan*  
Name

QC

Title

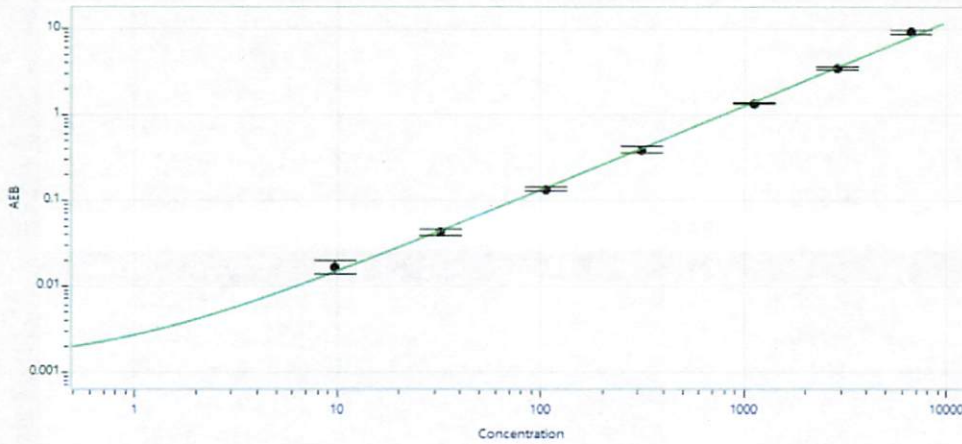
*Sonia Khillan 08SEP2023*  
Signature/Date

**NF-light® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	0.677
C	1.84
D	5.70
E	15.6
F	52.2
G	139
H	394

**GFAP® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	10.0
C	30.5
D	108
E	316
F	1115
G	3008
H	7350