

Product Number:	103520
Lot Number:	503409
Expiration:	20-Mar-2023
Platform(s):	HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103510	220610
Detector Reagent	103512	220611
SBG Reagent	103514	207903
Sample Diluent	103516	207904
Calibrators	103517	221303/224903
Control 1	103518	221505
Control 2	103519	221505
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).



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

Release Materials			
NF-light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	221505	4.27	3.10-5.46
Control 2	221505	431	338-507
Panel 1	207003	14.0	10.7-16.1
Panel 2	207003	6.77	4.77-7.68
Panel 3	207003	136	101-151
Panel 4	207003	388	273-416
GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	221505	86.3	72.2-112
Control 2	221505	10333	8148-12222
Panel 1	207003	60.1	49.4-83.9
Panel 2	207003	75.8	52.9-104
Panel 3	207003	910	684-1104
Panel 4	207003	6073	4259-7215

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

Review/Approval

*Sonia Khillan*

Name

*QC Manager*

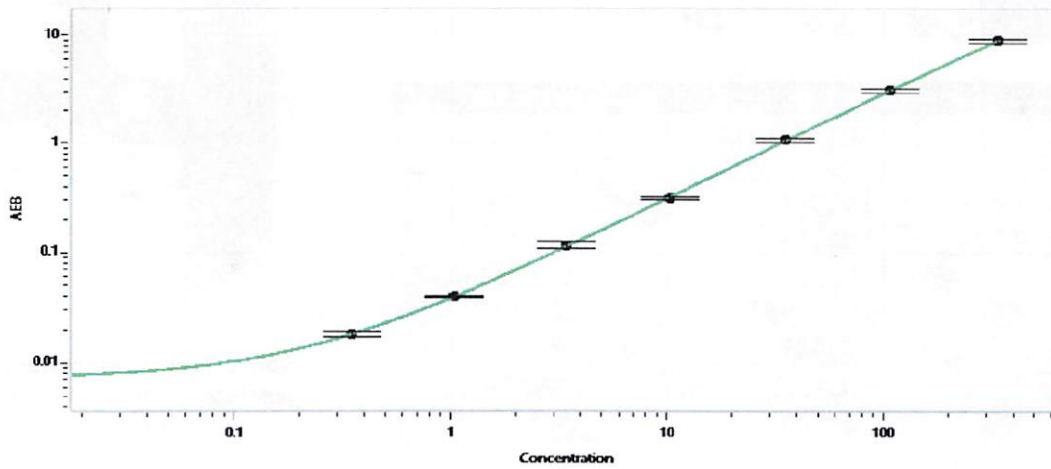
Title

*Sonia Khillan 1/16 SEP 2022*

Signature/Date

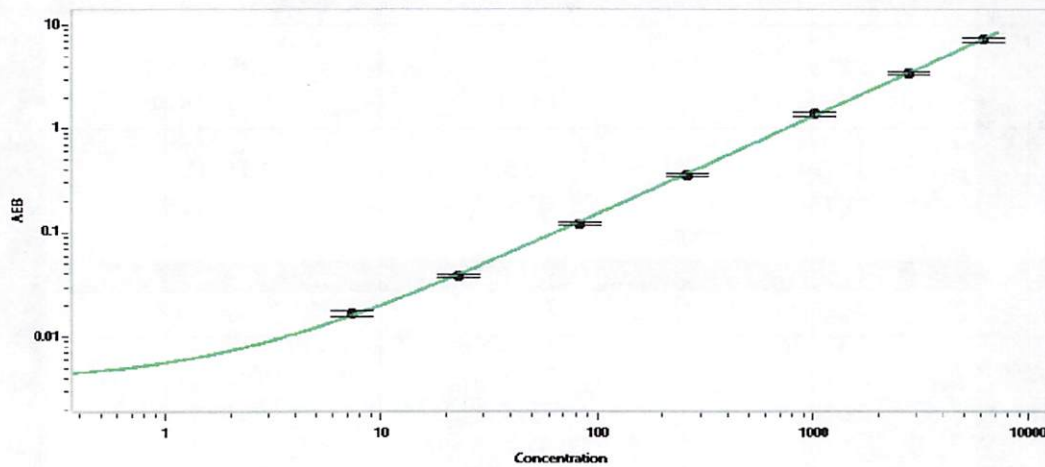


**NF-light® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	0.352
C	1.04
D	3.42
E	10.3
F	35.3
G	107
H	335

**GFAP® Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	7.36
C	22.9
D	82.5
E	261
F	1014
G	2766
H	6154