

Product Number:	103520
Lot Number:	503727
Expiration:	7-Sep-2023
Platform(s):	HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103510	227108
Detector Reagent	103512	227224
SBG Reagent	103514	227225
Sample Diluent	103516	227226
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>



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

<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	221505	4.07	2.87-5.47
Control 2	221505	437	335-502
Panel 1	207003	13.4	10.7-16.1
Panel 2	207003	6.96	4.77-7.68
Panel 3	207003	138	101-151
Panel 4	207003	389	273-416

**GFAP®**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	221505	97.2	73.6-111
Control 2	221505	10493	8061-12092
Panel 1	207003	59.2	49.4-83.9
Panel 2	207003	69.8	52.9-104
Panel 3	207003	805	684-1104
Panel 4	207003	5322	4259-7215

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

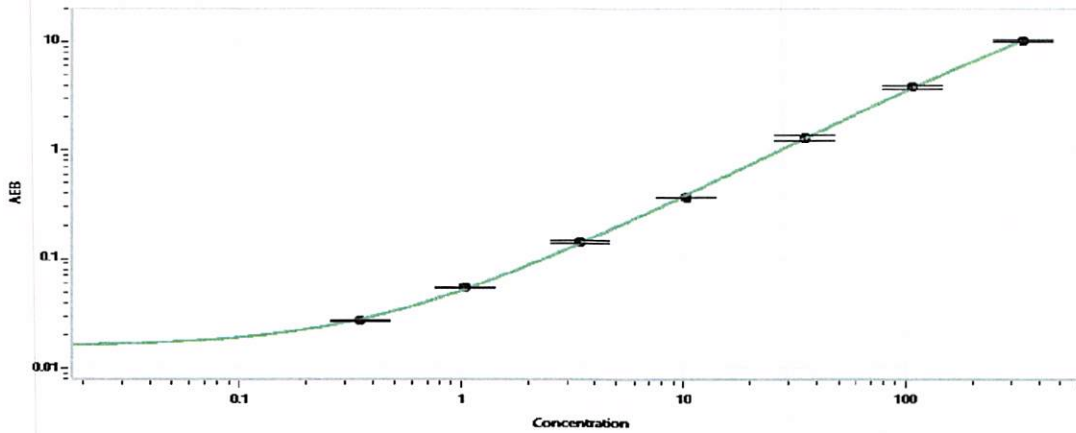
Review/Approval

*Sonia Khilari*  
Name

*QC Manager*  
Title

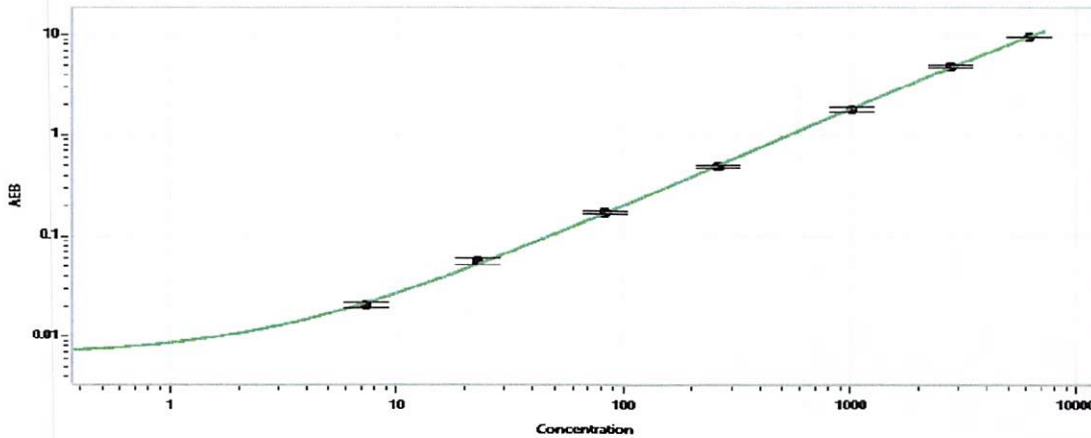
*Sonia Khilari - 10/08/2022*  
Signature/Date

**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