

Product Number:	103520
Lot Number:	504071
Expiration:	2-Sep-2024
Platform(s):	SR-X and HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103510	333510
Detector Reagent	103512	333506
SBG Reagent	103514	330402
Sample Diluent	103516	328512
Calibrators	103517	321906
Control 1	103518	321907
Control 2	103519	321907
SBG Diluent (SRX)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific

† Specific to use on the SRX platform - refer to Neuro 2-Plex B SR-X Kit Instructions (KI-0062).

Release Materials*			
NF-Light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	321907	5.55	4.78 - 7.17
Control 2	321907	553	439 - 658
Panel 1	310007	6.13	4.33 - 10.6
Panel 2	323304	5.24	4.46 - 8.31
Panel 3	306904	79.2	74.0 - 111
Panel 4	306904	462	359 - 597

GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	321907	82.3	60.0 - 116
Control 2	321907	8855	6348 - 10657
Panel 1	310007	65.3	53.5 - 103
Panel 2	323304	49.3	34.3 - 82.1
Panel 3	306904	1211	795 - 1440
Panel 4	306904	8661	6199 - 9299

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

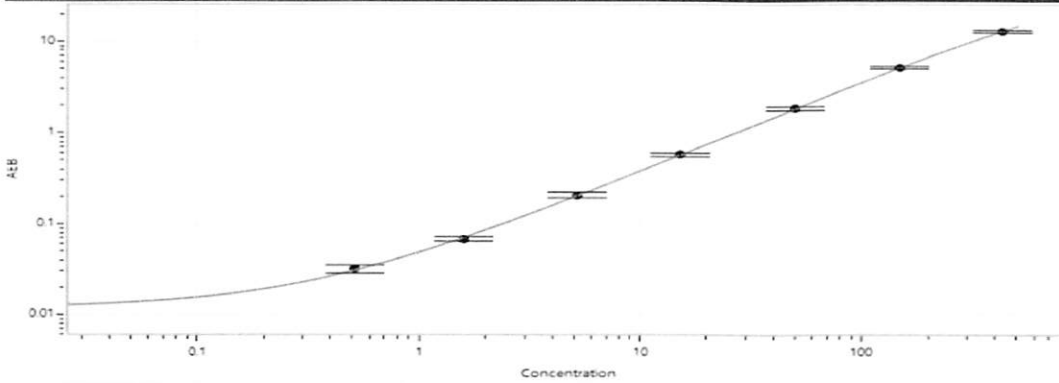
Review/Approval

*Sonia Khillan*  
Name

*Senior QC Analyst*  
Title

*Sonia Khillan 12-DEC-2023*  
Signature/Date

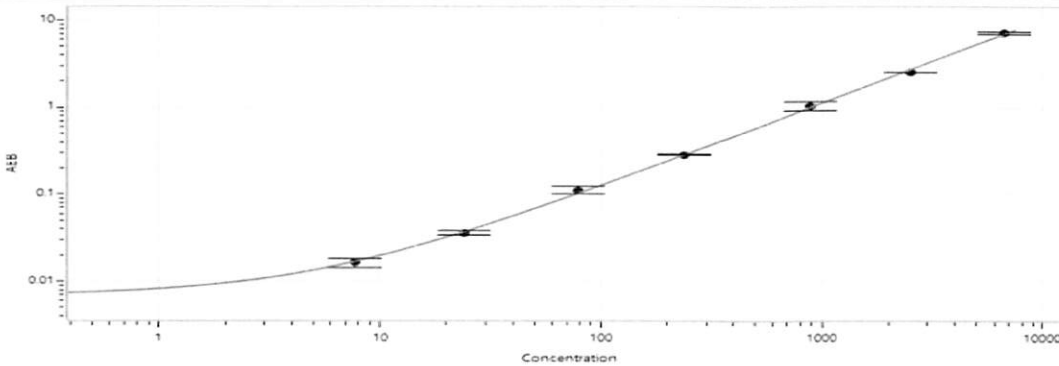
**NF-Light® Calibrator Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.517
C	1.59
D	5.17
E	15.1
F	50.4
G	149
H	431

**GFAP® Calibrator Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	7.66
C	24.2
D	78.8
E	237
F	880
G	2507
H	6673