

Product Number:	103520
Lot Number:	504042
Expiration:	5-Dec-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	333903
Detector Reagent	103512	334004
SBG Reagent	103514	334607
Sample Diluent	103516	334005
Calibrators	103517	321906
Control 1	103518	321907
Control 2	103519	321907
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ 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	5.67	4.72 - 7.13
Control 2	321907	530	436 - 655
Panel 1	310007	6.05	4.33 - 10.6
Panel 2	323304	5.20	4.46 - 8.31
Panel 3	306904	78.5	74.0 - 111
Panel 4	306904	434	359 - 597

GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	321907	89.0	61.6 - 114
Control 2	321907	8924	6504 - 10641
Panel 1	310007	76.9	53.5 - 103
Panel 2	323304	62.9	34.3 - 82.1
Panel 3	306904	1238	795 - 1440
Panel 4	306904	8515	6199 - 9299

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

Review/Approval

ROBERT SOMOGYI

QC MANAGER

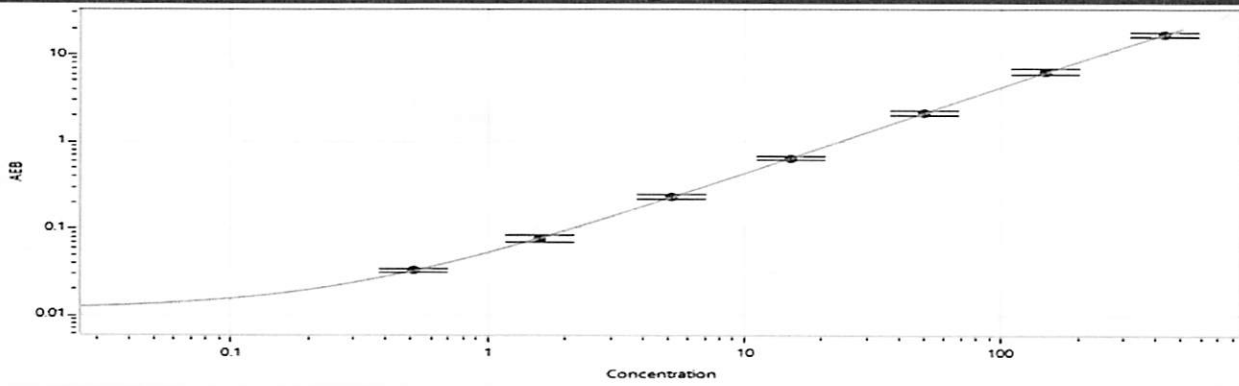
[Signature] 20 DEC 2023

Name

Title

Signature/Date

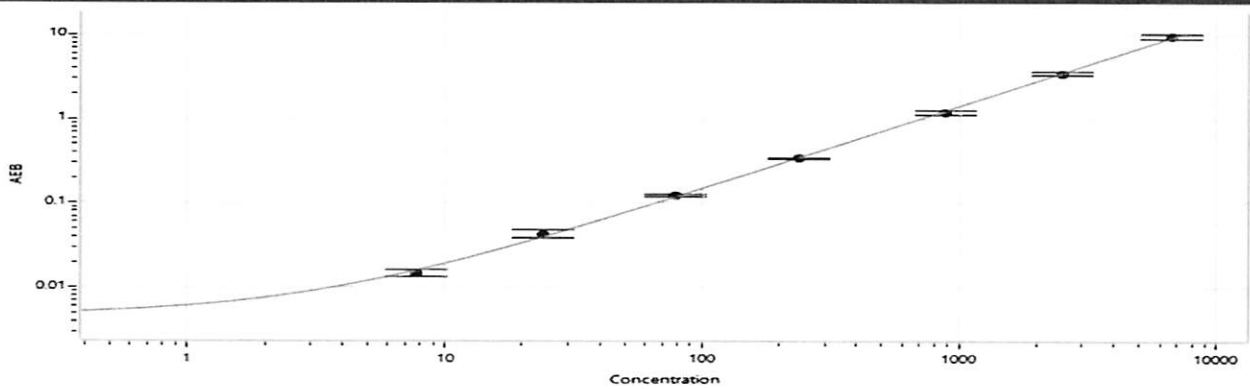
NF-Light® Calibrator Curve



**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