

Product Number:	103520
Lot Number:	504131
Expiration:	5-Aug-2025
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	403722
Detector Reagent	103512	403809
SBG Reagent	103514	408009
Sample Diluent	103516	403811
Calibrators	103517	403714
Control 1	103518	403715
Control 2	103519	403715
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	403715	5.79	4.06 - 6.67
Control 2	403715	472	384 - 577
Panel 1	405908	19.5	15.0 - 25.5
Panel 2	323304	6.36	4.46 - 8.31
Panel 3	306904	92.8	74.0 - 111
Panel 4	306904	469	359 - 597

GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	403715	87.8	68.8 - 111
Control 2	403715	9088	7370 - 11054
Panel 1	405908	119	74.8 - 150
Panel 2	323304	51.6	34.3 - 82.1
Panel 3	306904	1056	795 - 1440
Panel 4	306904	6822	6199 - 9299

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

Review/Approval

Cody Carlson

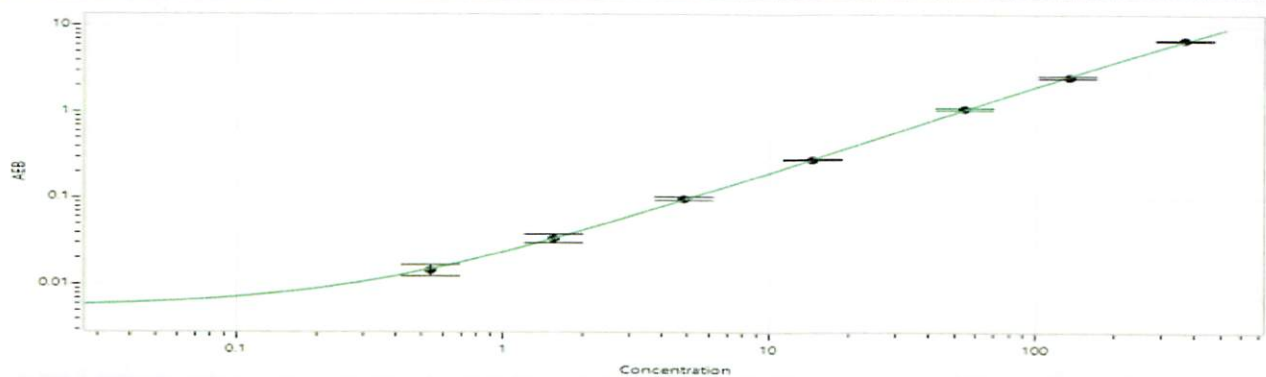
QC Manager

Name

Title

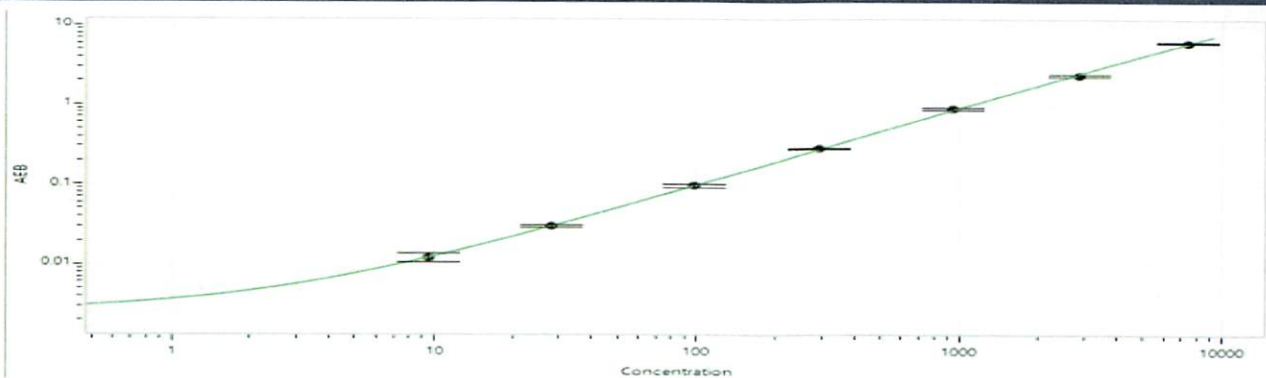
Cody Carlson 22 MAR 2024
Signature/Date

NF-Light® Calibrator Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.536
C	1.55
D	4.84
E	14.6
F	54.3
G	134
H	369

GFAP® Calibrator Curve



Calibrator Levels (pg/mL)	
A	0.000
B	9.46
C	27.9
D	97.3
E	292
F	937
G	2851
H	7379