

Product Number:	103520
Lot Number:	504137
Expiration:	3-Oct-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	409402
Detector Reagent	103512	409501
SBG Reagent	103514	409405
Sample Diluent	103516	409503
Calibrators	103517	413102
Control 1	103518	413103
Control 2	103519	413103
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	413103	5.13	4.09 - 6.57
Control 2	413103	585	468 - 702
Panel 1	405908	15.7	12.3 - 25.5
Panel 2	323304	5.27	4.46 - 8.31
Panel 3	410001	144	126 - 189
Panel 4	412411	695	543 - 814

GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	413103	93.0	67.8 - 120
Control 2	413103	9307	7344 - 11016
Panel 1	405908	133	76.1 - 165
Panel 2	323304	49.9	34.3 - 82.1
Panel 3	410001	467	346 - 519
Panel 4	412411	4422	3630 - 5445

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

Review/Approval

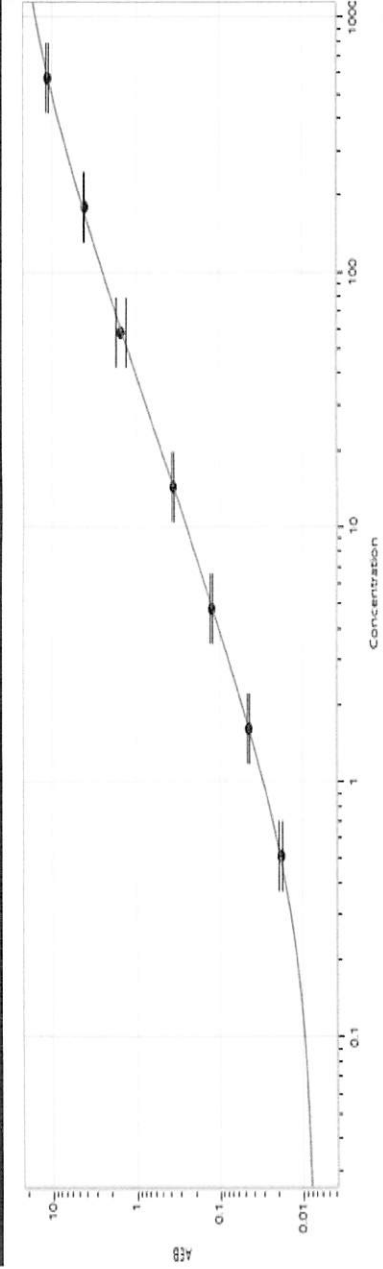
Cody Carlson

QC Manager

Name

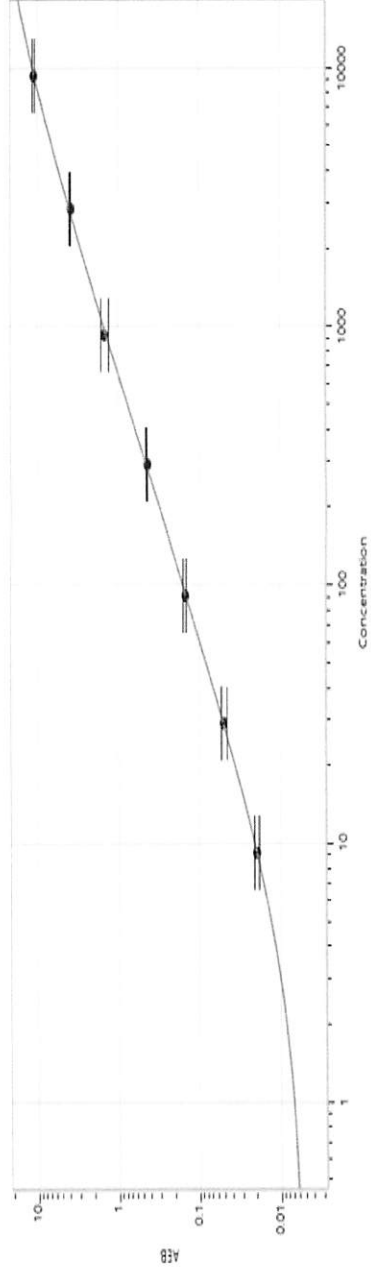
Title

[Signature] 11 JUN 2024
Signature/Date

NF-Light® Calibrator Curve

**Calibrator Levels
(pg/mL)**

A 0.000
 B 0.511
 C 1.62
 D 4.78
 E 14.4
 F 58.2
 G 178
 H 578

GFAP® Calibrator Curve

**Calibrator Levels
(pg/mL)**

A 0.000
 B 9.20
 C 29.0
 D 91.1
 E 294
 F 920
 G 2834
 H 9268