

Product Number:	103345
Lot Number:	504005
Expiration:	3-Nov-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	103338	330706
Detector Reagent	103339	331009
SBG Reagent	103340	335207
Sample Diluent	103341	331101
Calibrators	103342	330001
Control 1	103343	330002
Control 2	103344	330002
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

† Specific to use on the SR-X platform - refer to Neuro 4-Plex B Kit Instruction SR-X (KI-0123).

Review/Approval

Cody Carlson

QC Manager

[Signature] 21 Dec 2023

Name

Title

Signature/Date

Release Materials*

Tau

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	330002	2.91	2.09 - 3.44
Control 2	330002	106	80.6 - 130
Panel 2	306108	1.92	1.31 - 4.10
Panel 3	306108	24.2	20.8 - 31.2
Panel 4	306108	92.2	75.5 - 113

NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	330002	8.05	5.79 - 10.0
Control 2	330002	504	435 - 653
Panel 2	306108	12.8	9.47 - 16.5
Panel 3	306108	66.2	61.0 - 94.0
Panel 4	306108	286	255 - 473

GFAP®

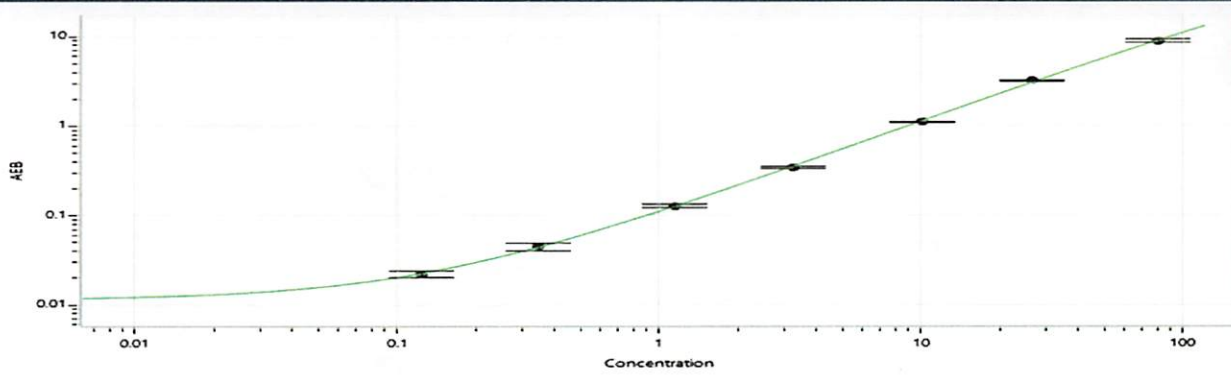
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	330002	289	215 - 323
Control 2	330002	12910	10183 - 15275
Panel 2	306108	102	52.2 - 127
Panel 3	306108	1154	751 - 1288
Panel 4	306108	6567	4770 - 7496

UCHL1®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	330002	159	123 - 184
Control 2	330002	13016	10165 - 15247
Panel 2	306108	119	62.2 - 195
Panel 3	306108	1777	1540 - 2321
Panel 4	306108	12061	10196 - 15294

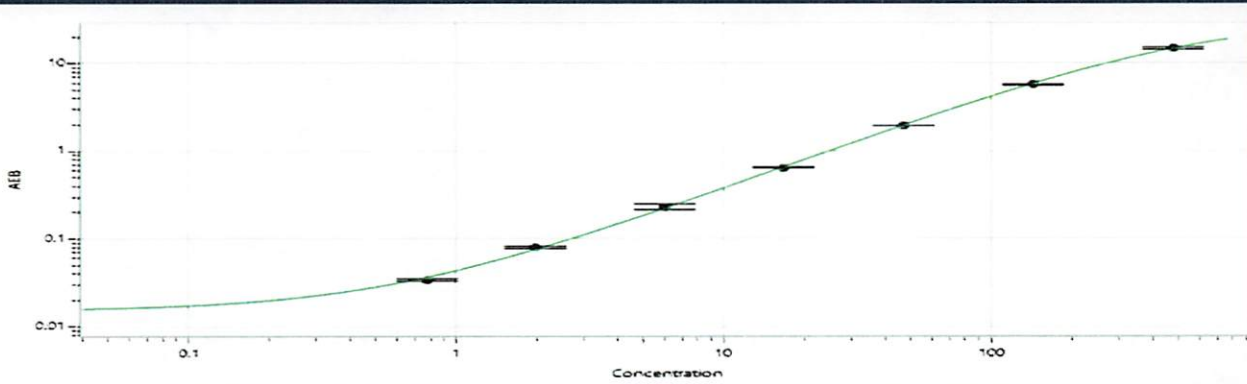
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Tau Calibrator Curve



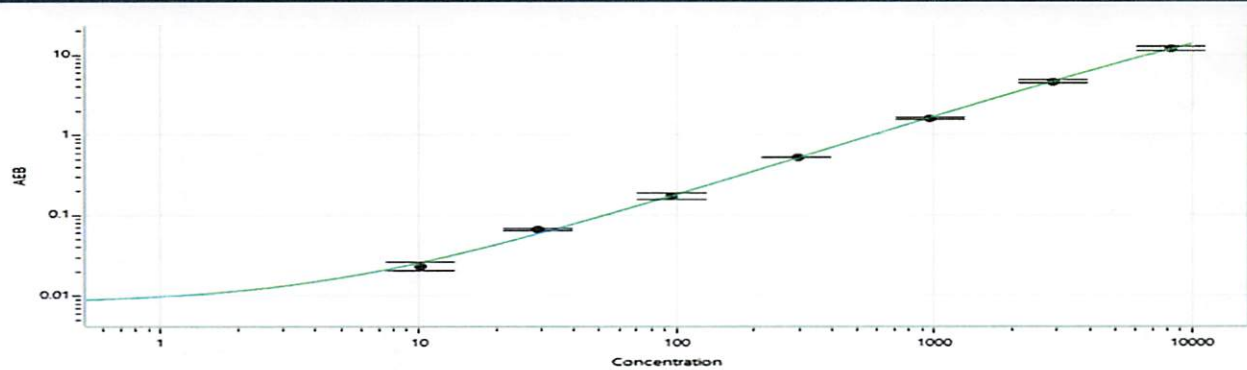
Calibrator Levels (pg/mL)	
A	0.000
B	0.124
C	0.347
D	1.15
E	3.23
F	10.1
G	26.7
H	80.1

NF-Light® Calibrator Curve



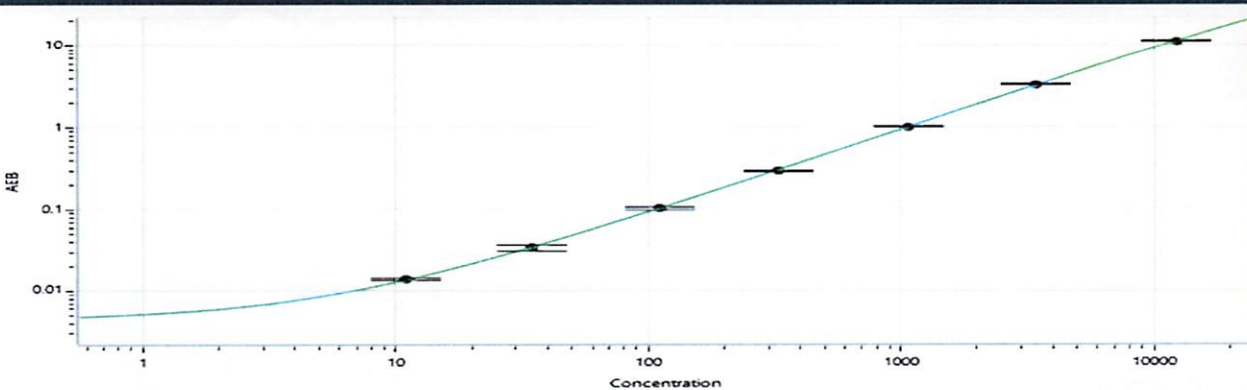
Calibrator Levels (pg/mL)	
A	0.000
B	0.784
C	1.97
D	6.04
E	16.6
F	46.7
G	143
H	475

GFAP® Calibrator Curve



Calibrator Levels (pg/mL)	
A	0.000
B	10.1
C	29.0
D	96.4
E	294
F	956
G	2890
H	8240

UCHL1® Calibrator Curve



Calibrator Levels (pg/mL)	
A	0.000
B	11.0
C	34.7
D	110
E	325
F	1060
G	3430
H	12310