

Product Number:	103345
Lot Number:	504312
Expiration:	3-May-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	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

Name

Title

Cody Carlson 12 Jul 2024
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

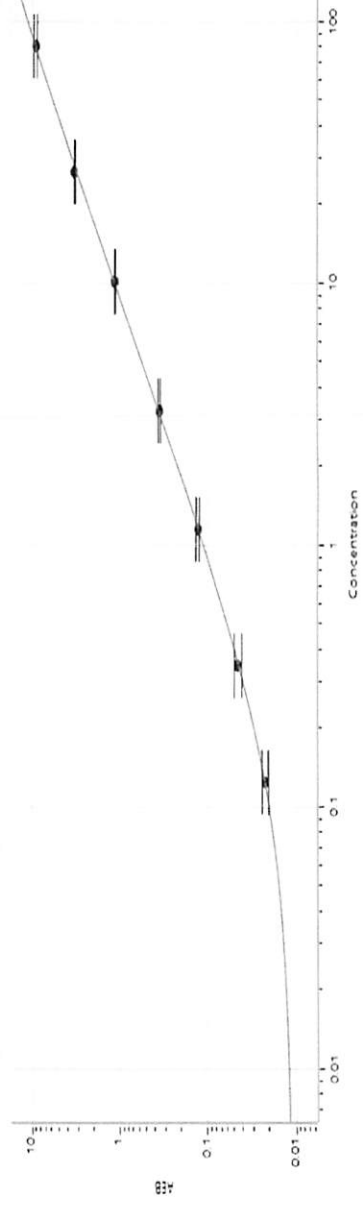
GFAP®

	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®

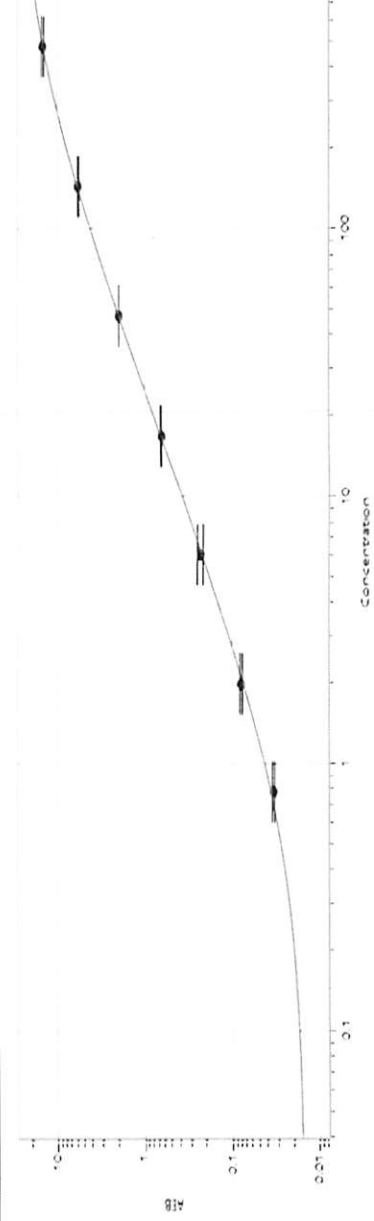
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	330002	159	123 - 184
Control 2	330002	13016	10165 - 15247
Panel 2	306108	119	62.2 - 195
Panel 3	306108	1777	1540 - 2321
Panel 4	306108	12061	10196 - 15294

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

Tau Calibrator Curve

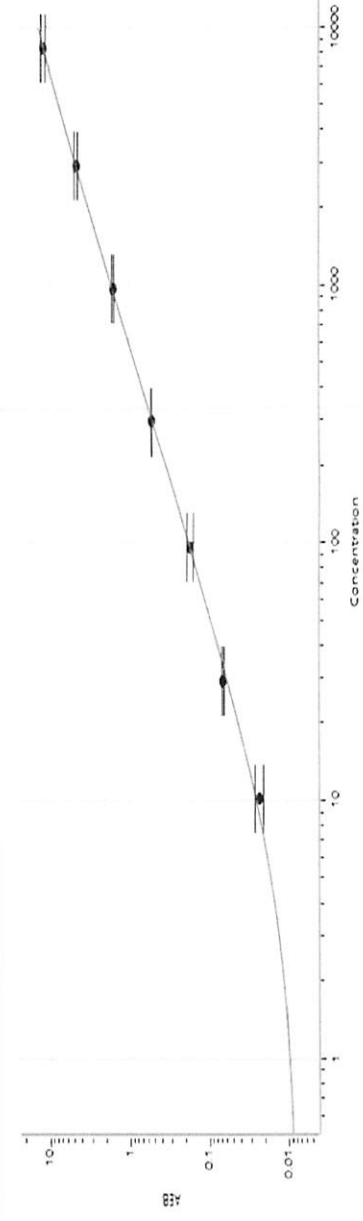
**Calibrator Levels
(pg/mL)**

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

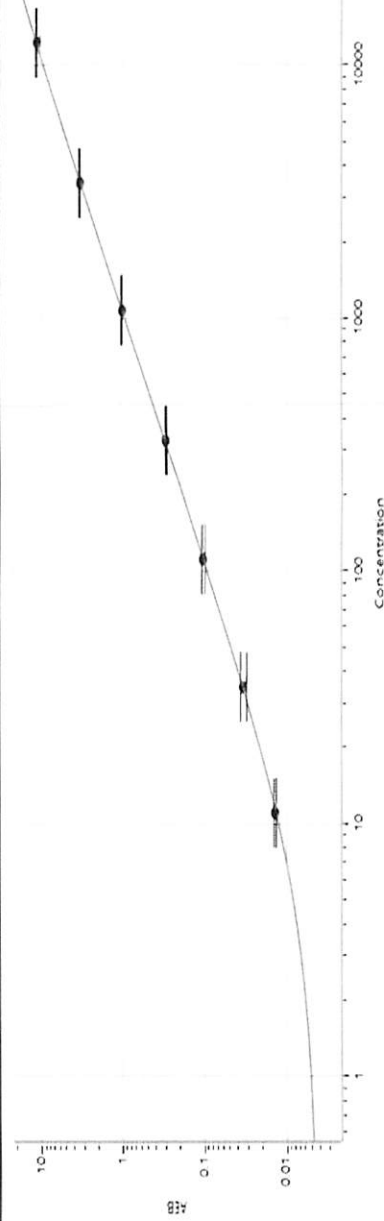
**Calibrator Levels
(pg/mL)**

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