

Product Number:	102153
Lot Number:	504167
Expiration:	2-Sep-2024
Platform(s):	SR-X and HD-X

Data below represents results generated on the Simoa® HD-X & SR-X Analyzer.

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	102141	325804	N/A
Detector Reagent	102143	325901	N/A
SBG Reagent	102145	328214	328214A
Sample Diluent	102147	326306	N/A
Calibrators	102150	325505	N/A
Control 1	102151	325506	N/A
Control 2	102152	325506	N/A
SBG Diluent (SR-X)†	103158	N/A ¹	N/A ¹
RGP Reagent	103159	N/A ¹	N/A ¹



¹ Reagents are not Kit Lot Specific

† Specific to use on the SRX platform - refer to Neuro 4-Plex A Kit Instruction SR-X (KI-0031).

Review/Approval

Cody Carlson

QC Manager

Name

Title

Cody Carlson 08MAR2024
Signature/Date

Release Materials*

Tau

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	325506	2.40	1.98 - 2.97
Control 2	325506	97.6	77.6 - 116
Panel 1	308106	31.6	25.3 - 37.9
Panel 2	309707	2.59	2.06 - 3.37
Panel 3	311701	83.4	59.2 - 88.8

NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	325506	5.66	4.54 - 8.09
Control 2	325506	558	470 - 705
Panel 1	308106	9.52	7.63 - 13.2
Panel 2	309707	7.39	6.02 - 9.14
Panel 3	311701	469	354 - 597

GFAP®

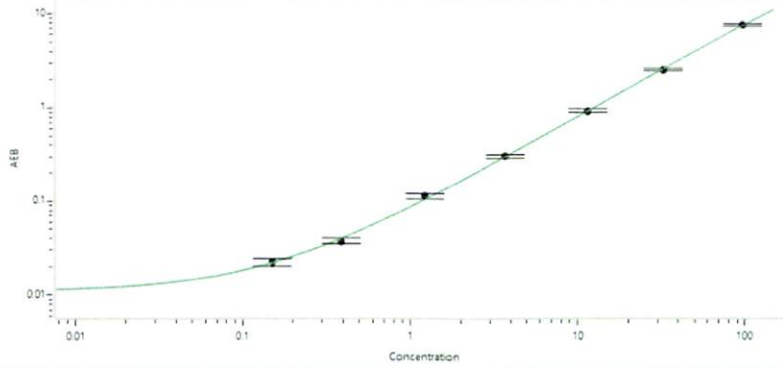
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	325506	26.0	20.0 - 30.1
Control 2	325506	1293	995 - 1492
Panel 1	308106	51.5	32.8 - 55.2
Panel 2	309707	90.7	64.9 - 97.4
Panel 3	311701	1119	872 - 1308

UCHL1®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	325506	144	88.8 - 181
Control 2	325506	11793	10006 - 15009
Panel 1	308106	94.4	40.6 - 142
Panel 2	309707	101	78.5 - 143
Panel 3	311701	5953	5347 - 8061

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

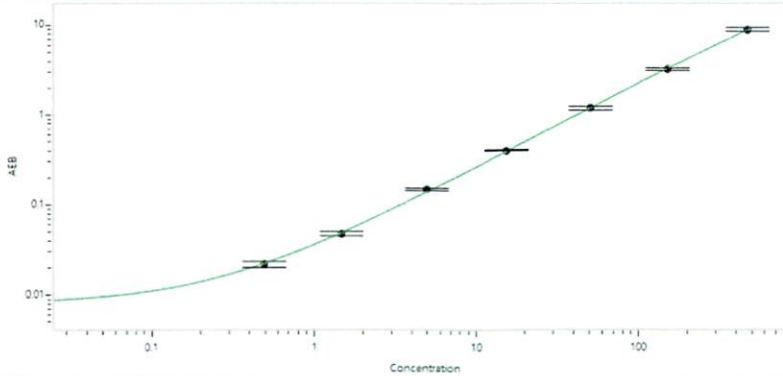
Tau Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.151
C	0.386
D	1.23
E	3.70
F	11.5
G	32.3
H	96.3

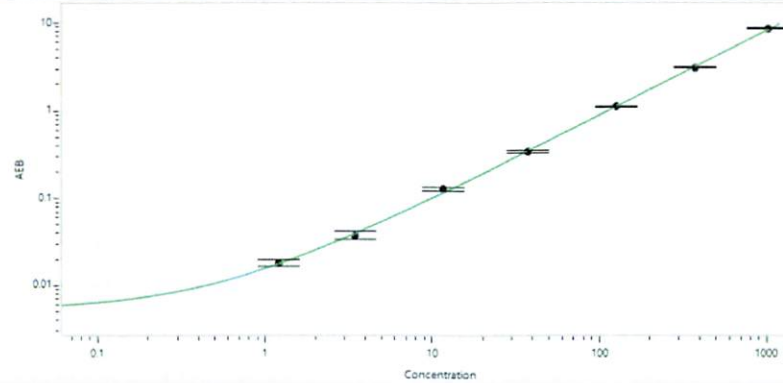
NF-Light® Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.496
C	1.48
D	5.00
E	15.5
F	50.4
G	153
H	467

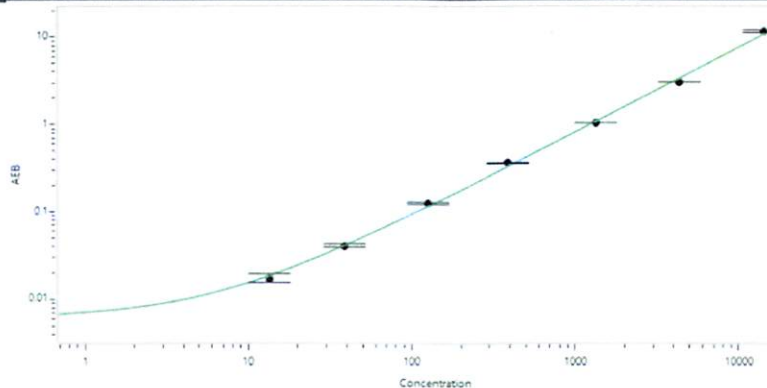
GFAP® Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	1.21
C	3.48
D	11.7
E	37.6
F	126
G	375
H	1019

UCHL1® Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	13.5
C	38.7
D	126
E	387
F	1337
G	4298
H	14194