

Product Number:	103345
Lot Number:	504050
Expiration:	09AUG2025
Platform(s):	SR-X and HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	103338	404006
Detector Reagent	103339	404306
SBG Reagent	103340	408001
Sample Diluent	103341	404307
Calibrators	103342	403902
Control 1	103343	403903
Control 2	103344	403903
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 SRX platform - refer to Neuro 4-Plex B Kit Instruction SR-X (KI-0123).



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

Review/Approval

Cody Carlson

QC Manager

Name

Title

Cody Carlson 12 APR 2024

Signature/Date

Release Materials*

Tau			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	403903	1.99	1.38 - 2.66
Control 2	403903	69.0	59.8 - 89.7
Panel 2	306108	2.02	1.31 - 4.06
Panel 3	306108	22.1	20.8 - 31.2
Panel 4	306108	83.4	75.5 - 113

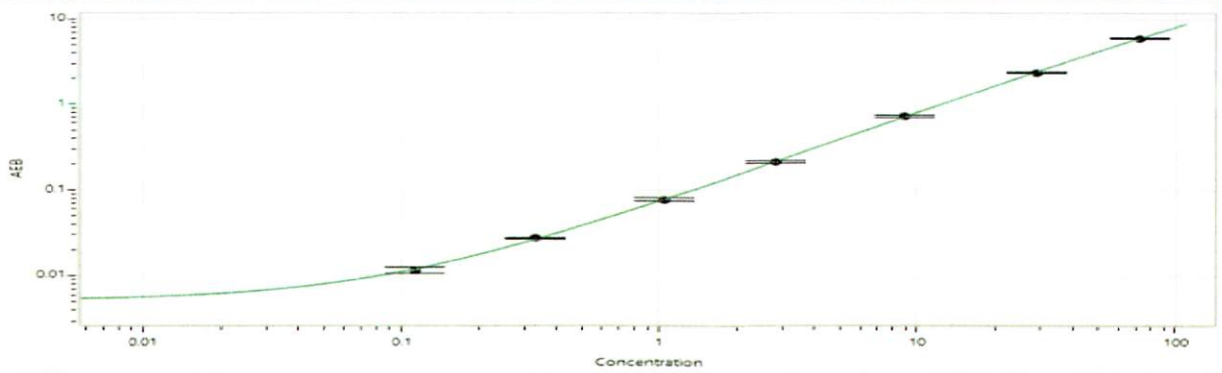
NF-Light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	403903	5.96	4.80 - 8.34
Control 2	403903	438	383 - 575
Panel 2	306108	11.9	9.47 - 16.5
Panel 3	306108	63.0	60.5 - 94.2
Panel 4	306108	276	255 - 473

GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	403903	97.2	77.8 - 135
Control 2	403903	9136	7695 - 11542
Panel 2	306108	85.1	52.2 - 127
Panel 3	306108	1099	751 - 1288
Panel 4	306108	6025	4770 - 7496

UCHL1®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	403903	135	108 - 161
Control 2	403903	11055	9165 - 13748
Panel 2	306108	120	62.2 - 195
Panel 3	306108	1759	1540 - 2321
Panel 4	306108	11690	10196 - 15294

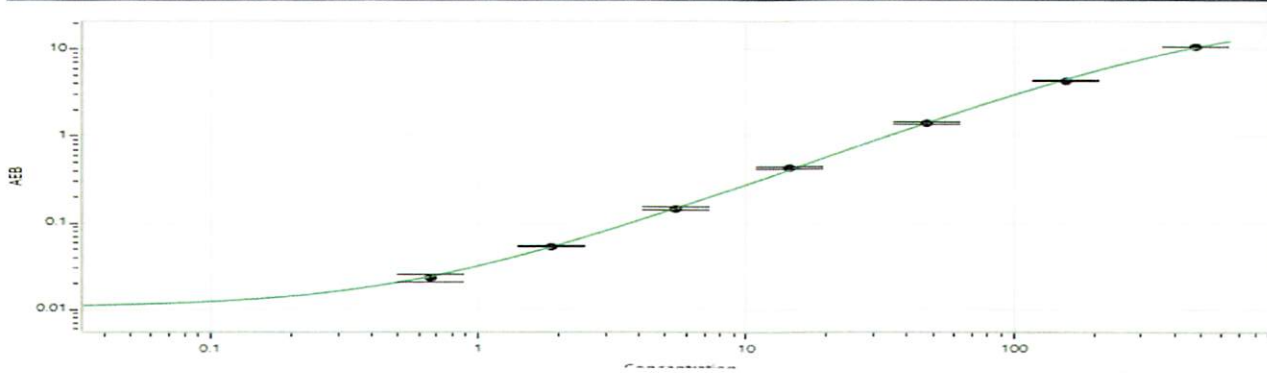
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Tau Calibrator Curve



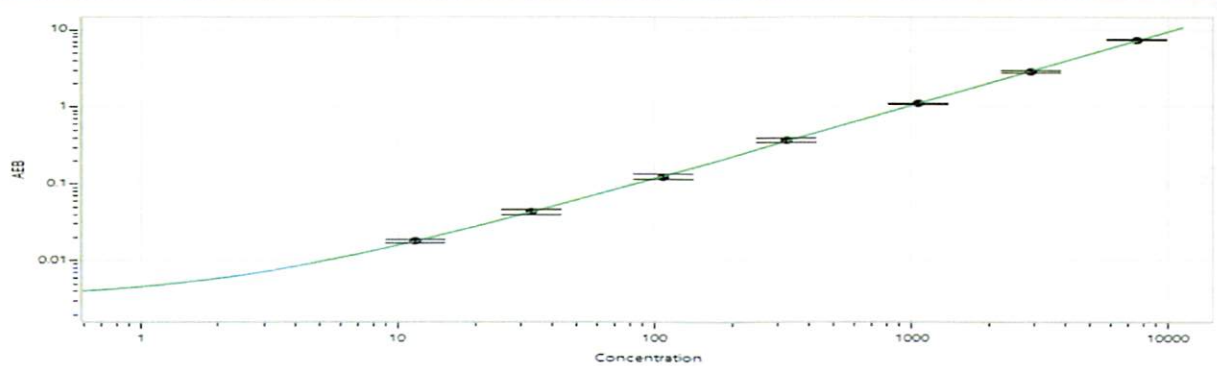
Calibrator Levels (pg/mL)	
A	0.000
B	0.113
C	0.330
D	1.04
E	2.82
F	8.87
G	29.2
H	72.6

NF-Light® Calibrator Curve



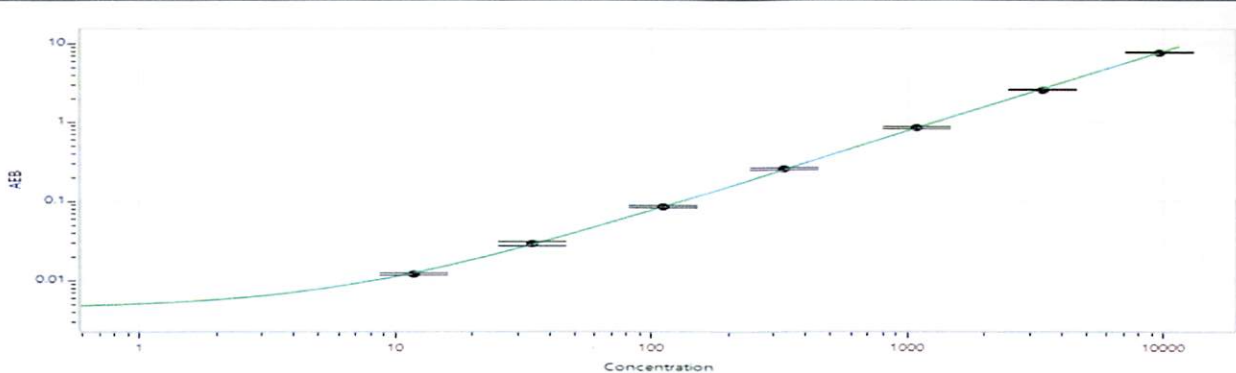
Calibrator Levels (pg/mL)	
A	0.000
B	0.661
C	1.87
D	5.45
E	14.6
F	47.4
G	157
H	482

GFAP® Calibrator Curve



Calibrator Levels (pg/mL)	
A	0.000
B	11.7
C	33.3
D	108
E	328
F	1072
G	2935
H	7545

UCHL1® Calibrator Curve



Calibrator Levels (pg/mL)	
A	0.000
B	11.7
C	34.1
D	112
E	329
F	1088
G	3349
H	9628