

Product Number:	102153
Lot Number:	504046
Expiration:	16-Aug-2025
Platform(s):	SR-X and HD-X

Component	Part Number	Lot Number
Bead Reagent	102141	404705
Detector Reagent	102143	417804
SBG Reagent	102145	408002
Sample Diluent Reagent	102147	405115
Reference Calibrators	102150	331901
Control 1	102151	331902
Control 2	102152	331902
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 A Kit Instruction SR-X (KI-0031).



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

Review/Approval

Cody Carlson

QC Manager

Name

Title

Cody Carlson 29 Jul 2024
Signature/Date

Release Materials*

Tau

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	331902	2.04	1.70 - 2.83
Control 2	331902	66.7	64.9 - 97.3

NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	331902	5.61	4.23 - 6.36
Control 2	331902	535	423 - 635

GFAP®

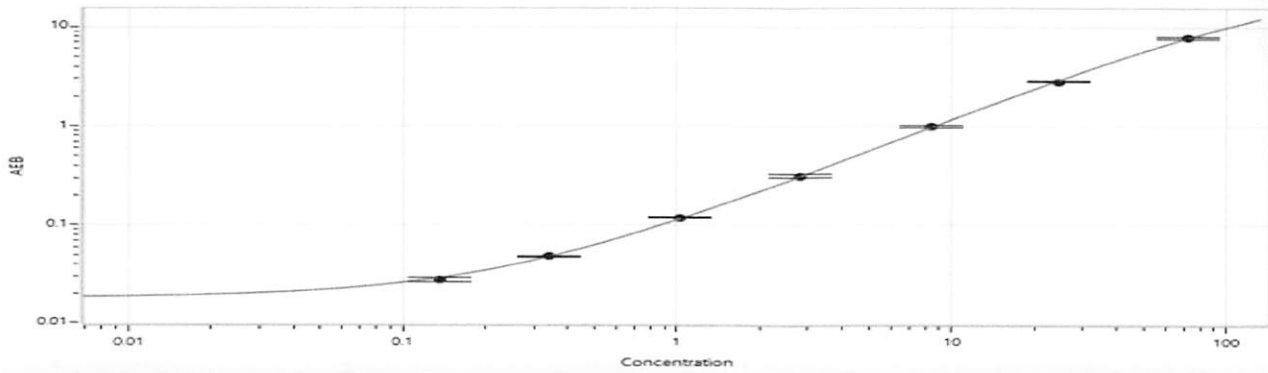
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	331902	20.2	14.7 - 22.1
Control 2	331902	1012	785 - 1178

UCHL1®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	331902	144	91.8 - 147
Control 2	331902	11006	8264 - 12397

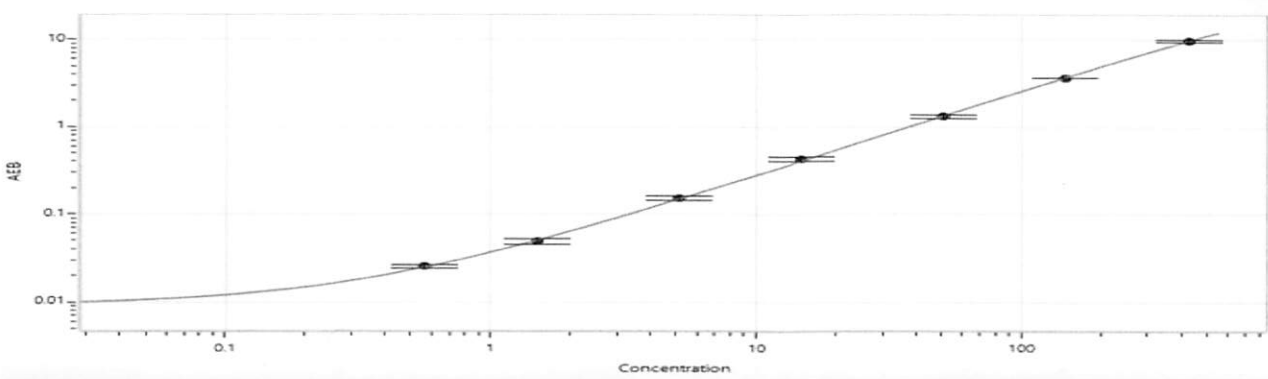
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Tau Calibrator Curve



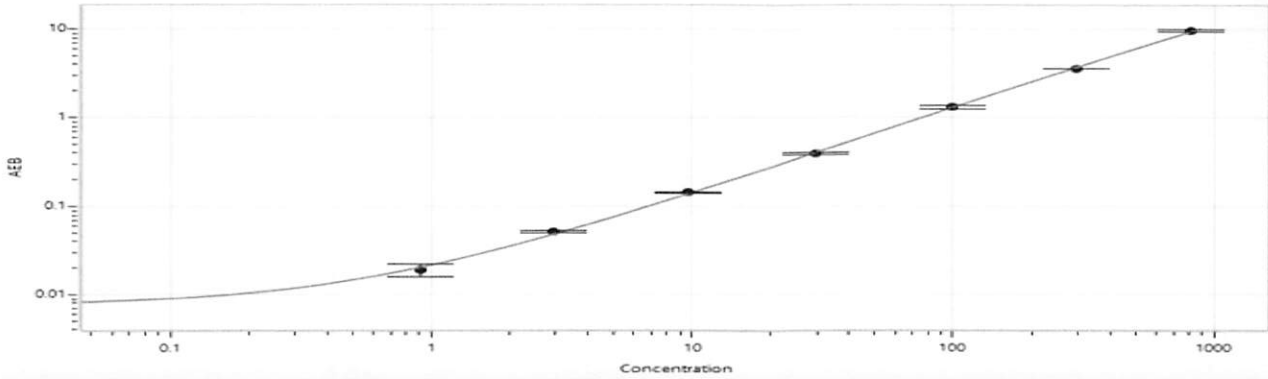
Calibrator Levels (pg/mL)	
A	0.000
B	0.136
C	0.339
D	1.02
E	2.79
F	8.45
G	24.4
H	72.5

NF-Light® Calibrator Curve



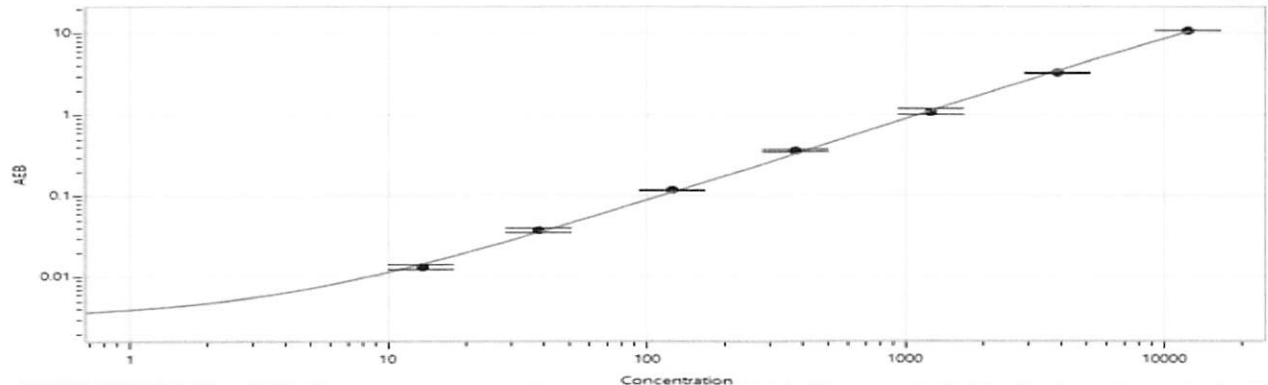
Calibrator Levels (pg/mL)	
A	0.000
B	0.565
C	1.51
D	5.16
E	14.9
F	51.0
G	147
H	426

GFAP® Calibrator Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.909
C	2.93
D	9.65
E	29.6
F	98.6
G	294
H	819

UCHL1® Calibrator Curve



Calibrator Levels (pg/mL)	
A	0.000
B	13.5
C	38.2
D	126
E	373
F	1257
G	3883
H	12464