

Product Number: 103345
Lot Number: 503475
Expiration: 18-Jun-2023
Platform(s): HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103338	216903
Detector Reagent	103339	220812
SBG Reagent	103340	217306
Sample Diluent	103341	217307
Calibrators	103342	225803
Control 1	103343	217802
Control 2	103344	217802
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

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



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

Release Materials

Tau			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	217802	2.10	1.65-2.63
Control 2	217802	78.9	64.7-106
Panel 1	206807	13.8	10.8-16.2
Panel 2	206807	9.26	6.79-10.2
Panel 3	206807	26.9	21.3-32.0
Panel 4	206807	90.2	74.7-112

NF-light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	217802	4.76	3.62-5.64
Control 2	217802	431	296-588
Panel 1	206807	9.62	7.65-11.5
Panel 2	206807	3.70	2.60-4.66
Panel 3	206807	131	101-161
Panel 4	206807	666	553-866

GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	217802	129	86.9-166
Control 2	217802	9503	6831-12274
Panel 1	206807	67.6	51.2-81.9
Panel 2	206807	29.6	11.9-36.0
Panel 3	206807	981	766-1149
Panel 4	206807	4804	3897-5845

UCH-L1®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	217802	107	81.8-135
Control 2	217802	9601	7924-11886
Panel 1	206807	89.6	37.5-102
Panel 2	206807	84.1	33.0-95.4
Panel 3	206807	1196	794-1250
Panel 4	206807	7570	5671-9023

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

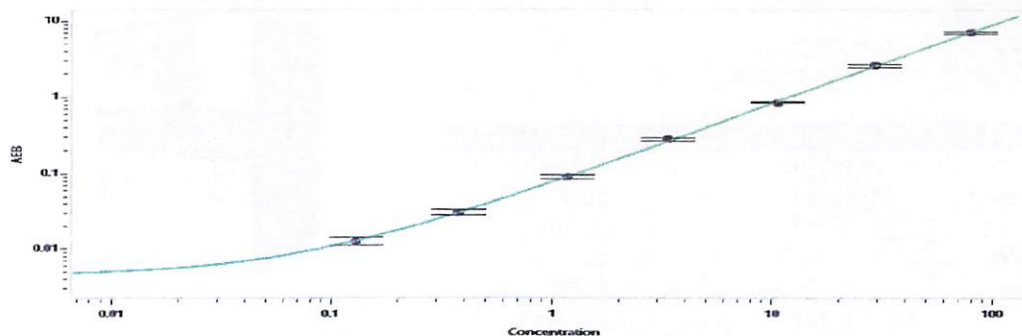
Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 30SEP2022
Signature/Date

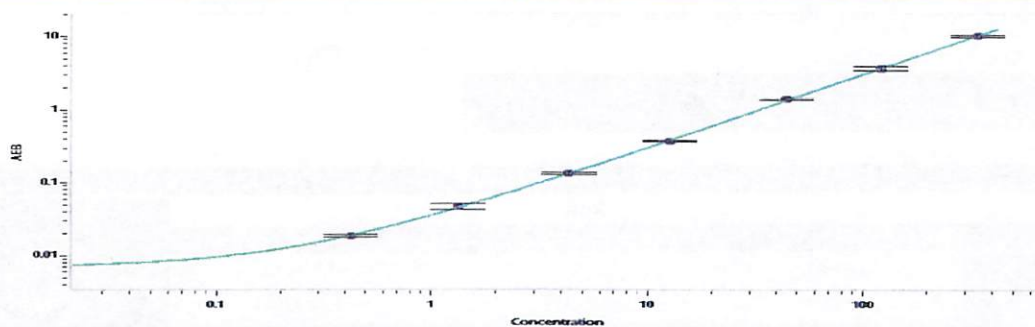
Tau Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.131
C	0.375
D	1.17
E	3.35
F	10.6
G	29.1
H	78.5

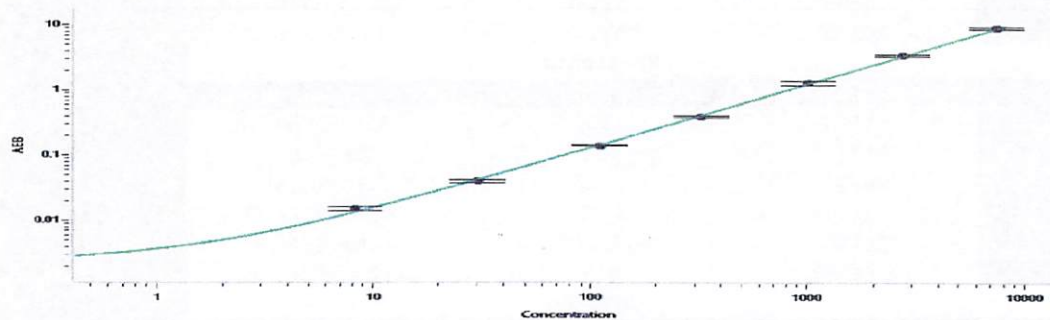
NF-light® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.427
C	1.330
D	4.36
E	12.7
F	44.9
G	122
H	340

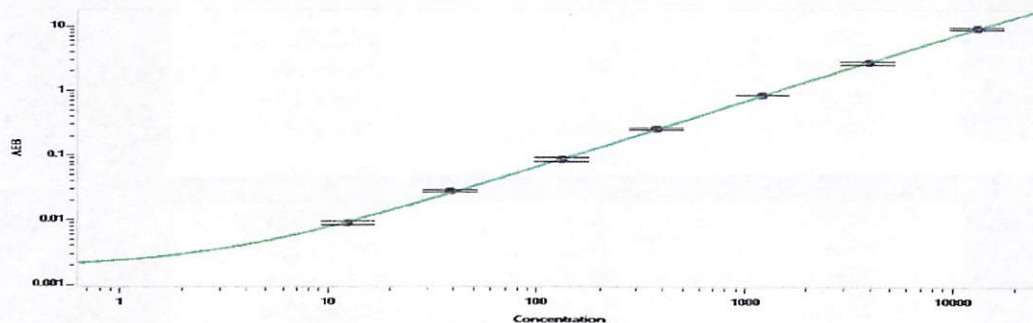
GFAP® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	8.22
C	30.3
D	109
E	320
F	1012
G	2741
H	7457

UCHL1® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	12.5
C	38.4
D	132
E	373
F	1223
G	3926
H	13117