

Product Number: 103345
Lot Number: 503202
Expiration: 16-Oct-2022
Platform(s): HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103338	130301
Detector Reagent	103339	131514
SBG Reagent	103340	130304
Sample Diluent	103341	130305
Calibrators	103342	131309
Control 1	103343	131310
Control 2	103344	131310
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)	Mean Range (pg/mL)
Control 1	131310	2.63	1.62 - 2.97
Control 2	131310	108	63.3 - 126
Panel 1	111709	5.12	4.40-7.25
Panel 2	111709	4.49	4.13 - 7.17
Panel 3	111709	35.3	25.9 - 39.7
Panel 4	111709	96.0	49.9 - 128

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	131310	7.46	5.15 - 8.37
Control 2	131310	551	433 - 651
Panel 1	111709	5.73	3.34-6.62
Panel 2	111709	11.5	6.78 - 12.3
Panel 3	111709	190	130 - 230
Panel 4	111709	676	351 - 831

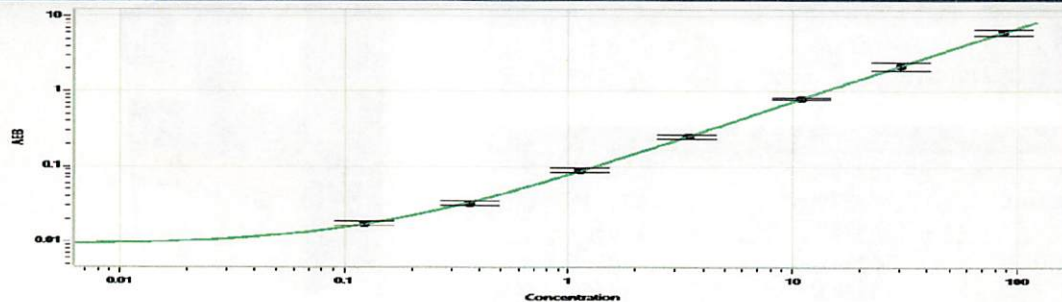
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	131310	176	115 - 206
Control 2	131310	10704	8431 - 12646
Panel 1	111709	36.3	21.5-54.5
Panel 2	111709	118	59.5 - 127
Panel 3	111709	1507	1268 - 1902
Panel 4	111709	10264	6727 - 13489

UCH-L1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	131310	163	118 - 182
Control 2	131310	11369	9537 - 14306
Panel 1	111709	94.2	62.0-129
Panel 2	111709	100	68.5 - 136
Panel 3	111709	1597	1352 - 2080
Panel 4	111709	6918	4654 - 9516

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

Review/Approval: Junpil Park, QC Supervisor
Name: Junpil Park Title: QC Supervisor Signature/Date: [Signature] 30 Nov 2021

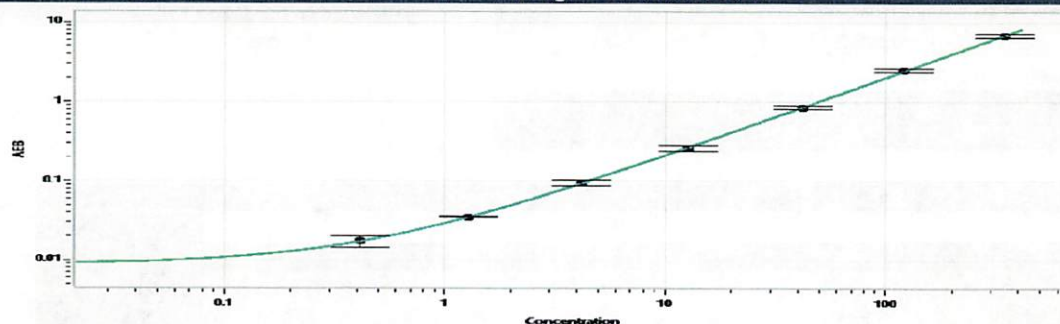
Tau Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.124
C	0.367
D	1.12
E	3.41
F	10.9
G	30.4
H	87.6

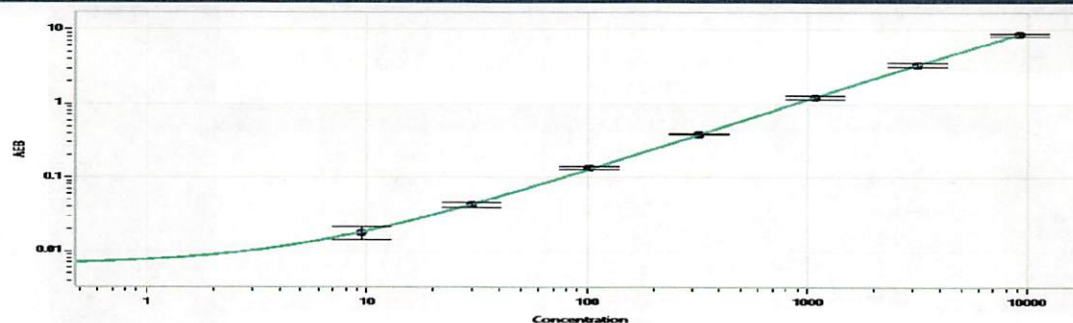
NF-light® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.418
C	1.28
D	4.18
E	12.8
F	42.0
G	120
H	346

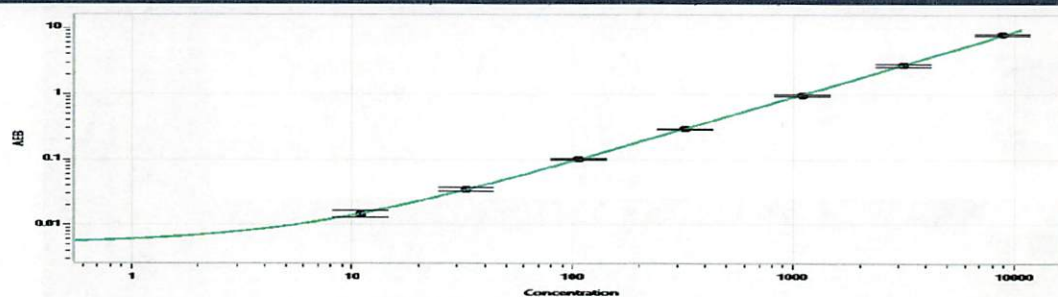
GFAP® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	9.42
C	29.9
D	102
E	323
F	1090
G	3175
H	9115

UCHL1® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	10.9
C	32.6
D	106
E	324
F	1104
G	3140
H	8825