

Product Number: 103345
Lot Number: 503228
Expiration: 3-Sep-2022
Platform(s): HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103338	132401
Detector Reagent	103339	133801
SBG Reagent	103340	132608
Sample Diluent	103341	132811
Calibrators	103342	131910
Control 1	103343	132304
Control 2	103344	132304
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	132304	1.95	1.48 - 2.34
Control 2	132304	83.6	53.4 - 94.6
Panel 1	111709	5.19	4.21 - 7.16
Panel 2	111709	6.32	4.20 - 7.40
Panel 3	111709	30.6	25.8 - 39.0
Panel 4	111709	91.1	54.8 - 124

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	132304	4.93	3.29 - 8.38
Control 2	132304	478	368 - 564
Panel 1	111709	4.26	3.12 - 6.52
Panel 2	111709	9.36	6.88 - 12.1
Panel 3	111709	188	134 - 229
Panel 4	111709	662	378 - 835

GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	132304	107	74.3 - 219
Control 2	132304	11908	8592 - 13132
Panel 1	111709	31.7	19.6 - 53.7
Panel 2	111709	88.0	58.5 - 125
Panel 3	111709	1830	1103 - 2104
Panel 4	111709	11899	6912 - 14100

UCH-L1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	132304	97.9	83.8 - 132
Control 2	132304	9172	7049 - 10574
Panel 1	111709	81.6	58.6 - 126
Panel 2	111709	107	73.1 - 134
Panel 3	111709	1767	1359 - 2096
Panel 4	111709	8384	4796 - 9951

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

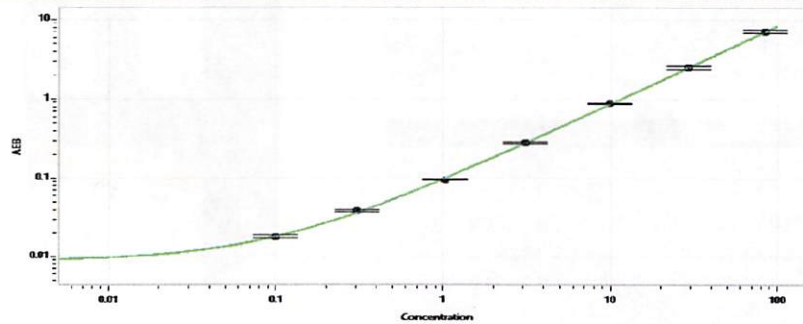
Review/Approval

Junpil Park
Name

QC Supervisor
Title

 15 Dec 2021
Signature/Date

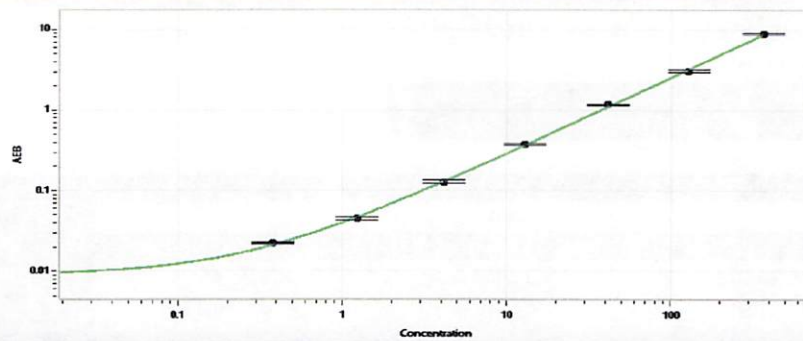
Tau Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.100
C	0.306
D	1.02
E	3.07
F	9.91
G	29.4
H	83.9

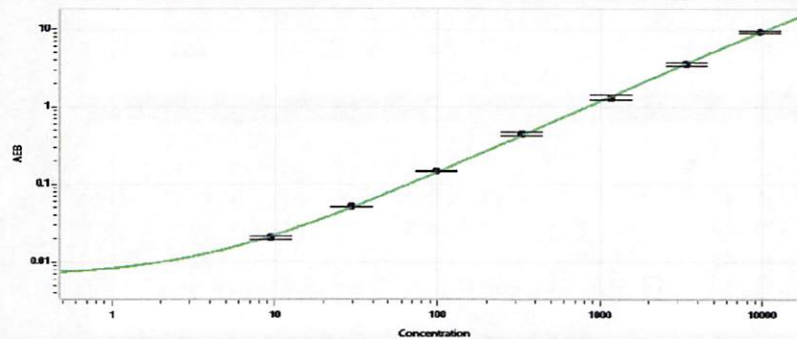
NF-light® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.376
C	1.23
D	4.13
E	12.9
F	41.7
G	128
H	366

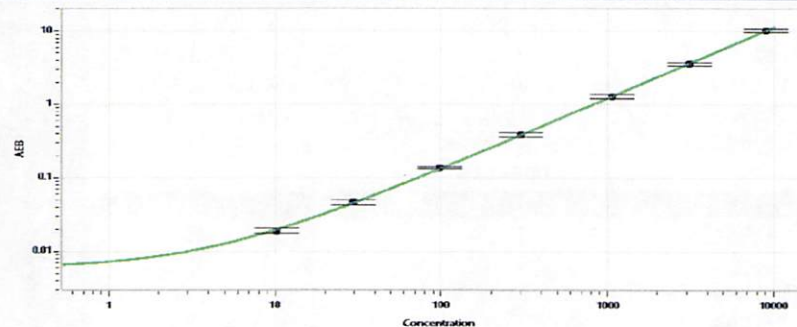
GFAP® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	9.41
C	29.5
D	98.2
E	325
F	1158
G	3388
H	9588

UCHL1® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	10.3
C	29.6
D	97.5
E	300
F	1060
G	3133
H	9008