

Product Number: 103345
Lot Number: 503266
Expiration: 3-Sep-2022
Platform(s): HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	103338	134707
Detector Reagent	103339	134706
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).

Release Materials

Tau			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	132304	1.79	1.47 - 2.29
Control 2	132304	76.0	55.8 - 93.0
Panel 1	111709	5.35	4.25 - 7.00
Panel 2	111709	6.27	4.37 - 7.41
Panel 3	111709	27.1	23.4 - 39.3
Panel 4	111709	84.0	56.7 - 120

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	132304	5.25	3.34 - 8.04
Control 2	132304	443	365 - 555
Panel 1	111709	5.00	3.21 - 6.49
Panel 2	111709	8.88	6.87 - 11.9
Panel 3	111709	178	135 - 227
Panel 4	111709	568	389 - 810

GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	132304	116	67.0 - 209
Control 2	132304	11059	8729 - 13094
Panel 1	111709	32.7	19.7 - 52.1
Panel 2	111709	92.2	60.7 - 123
Panel 3	111709	1587	1137 - 2064
Panel 4	111709	9848	7055 - 13718

UCH-L1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	132304	106	85.9 - 129
Control 2	132304	8854	7058 - 10587
Panel 1	111709	96.9	61.8 - 124
Panel 2	111709	112	75.7 - 135
Panel 3	111709	1745	1380 - 2082
Panel 4	111709	7120	4981 - 9673

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

Review/Approval

Junpil Park

Name

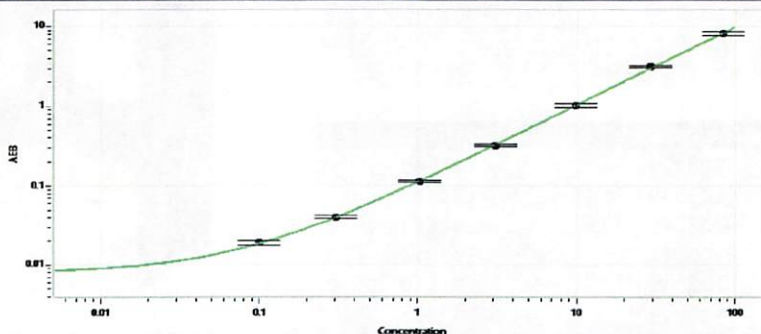
QC Supervisor

Title

Signature/Date

 20 Dec 2021

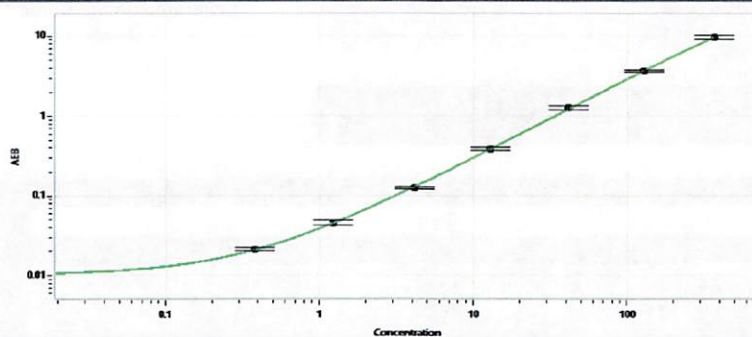
Tau Calibration Curve



**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

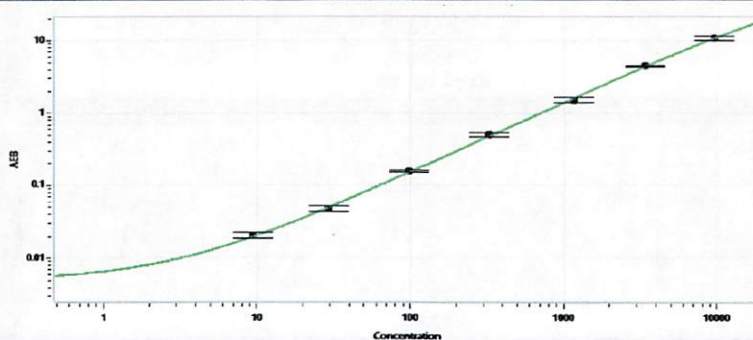
NF-light® Calibration Curve



**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

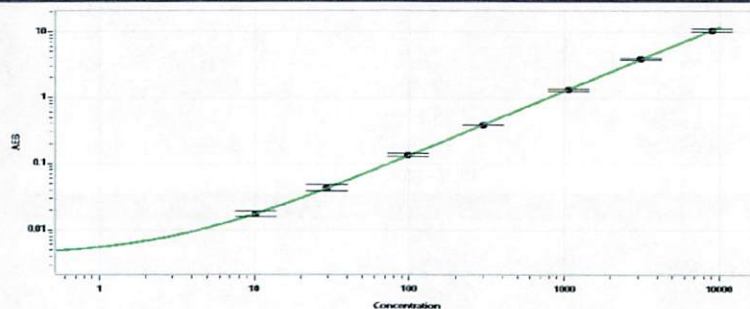
GFAP® Calibration Curve



**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