

Product Number:	103345
Lot Number:	502572
Expiration:	7-Oct-2021
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103338	028107
Detector Reagent	103339	028203
SBG Reagent	103340	028307
Sample Diluent	103341	028312
Calibrators	103342	033705
Control 1	103343	028901
Control 2	103344	028901
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	028901	1.74	1.22-1.84
Control 2	028901	90.8	63.7-107
Panel 2	921102	1.12	0.312 - 2.93
Panel 3	921103	26.0	20.3 - 34.0
Panel 4	921104	41.5	29.9 - 61.1

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	028901	8.26	6.75-11.1
Control 2	028901	932	697-1045
Panel 2	921102	6.98	5.75 - 8.63
Panel 3	921103	193	147 - 221
Panel 4	921104	591	555 - 832

GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	028901	111	77.6-116
Control 2	028901	12077	9096-13645
Panel 2	921102	42.9	27.8 - 53.1
Panel 3	921103	979	638 - 1195
Panel 4	921104	5337	3584 - 6789

UCH-L1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	028901	130	94.8-142
Control 2	028901	12167	8724-13086
Panel 2	921102	101	84.5 - 127
Panel 3	921103	3017	2398 - 4156
Panel 4	921104	7430	6436 - 10728

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

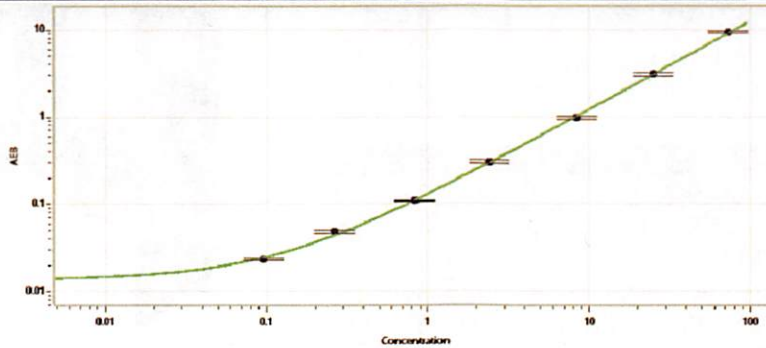
Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 14Dec 2020
Signature/Date

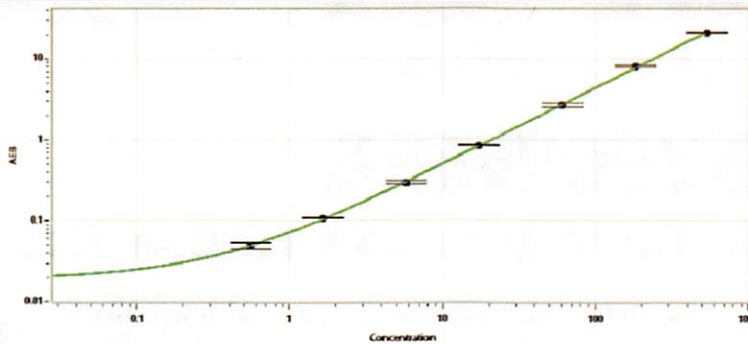
Tau Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.096
C	0.263
D	0.833
E	2.41
F	8.42
G	24.9
H	72.6

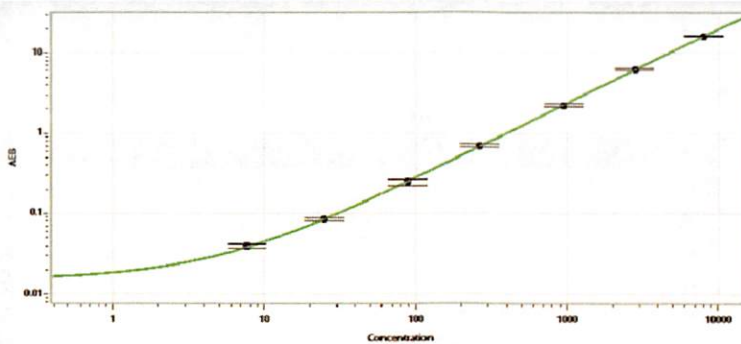
NF-light® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.552
C	1.64
D	5.78
E	17.2
F	60.7
G	181
H	535

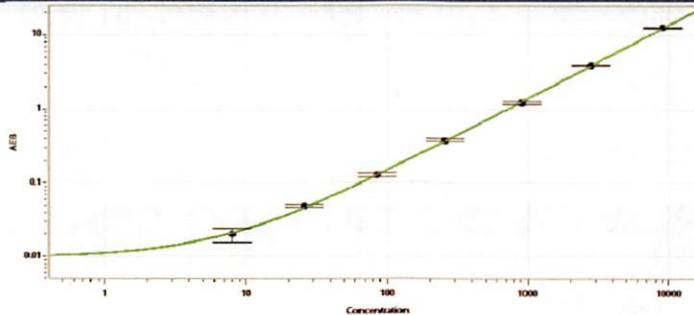
GFAP® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	7.72
C	24.9
D	87.8
E	261
F	951
G	2808
H	8016

UCHL1® Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	8.04
C	25.9
D	85.4
E	256
F	910
G	2778
H	9075