

Product Number:	103345
Lot Number:	502280
Expiration:	15-Jan-2021
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103338	006410
Detector Reagent	103339	006413
SBG Reagent	103340	006502
Sample Diluent	103341	006503
Calibrators	103342	931901
Controls	103343-103344	931902
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	931902	2.08	1.41-2.42
Control 2	931902	96.0	72.1-108
Panel 2	921102	0.912	0.557-2.95
Panel 3	921103	27.6	22.4-33.6
Panel 4	921104	45.9	36.4-58.8

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931902	6.18	4.86-7.29
Control 2	931902	543	439-658
Panel 2	921102	8.02	5.74-8.61
Panel 3	921103	205	145-217
Panel 4	921104	789	542-813

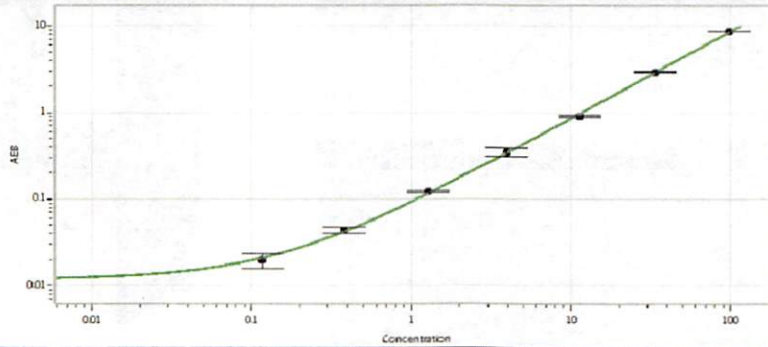
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931902	94.8	73.7-111
Control 2	931902	10531	8163-12244
Panel 2	921102	43.8	27.5-53.5
Panel 3	921103	997	625-1222
Panel 4	921104	6146	3418-6873

UCH-L1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931902	149	70.8-168
Control 2	931902	11911	8161-13158
Panel 2	921102	114	84.3-127
Panel 3	921103	3577	2320-4225
Panel 4	921104	9285	6239-10760

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

Review/Approval Patrizia Stadler Assoc. Director, QC p.stadler 23/11/2020
Name Title Signature/Date

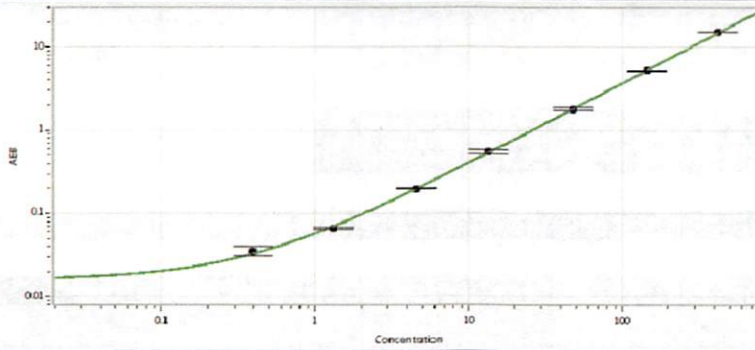
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.117
C	0.377
D	1.28
E	3.95
F	11.4
G	33.7
H	97.8

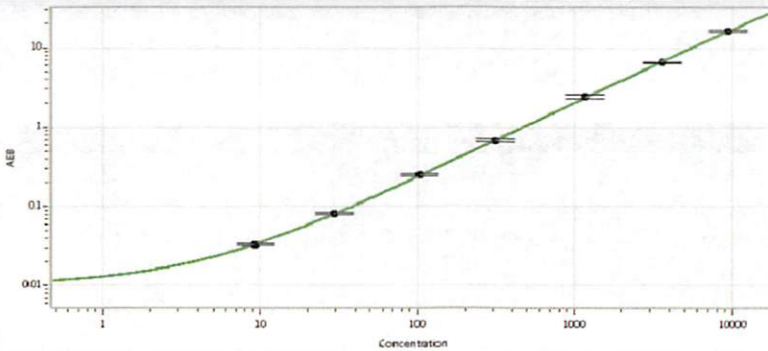
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.393
C	1.32
D	4.58
E	13.3
F	47.9
G	146
H	421

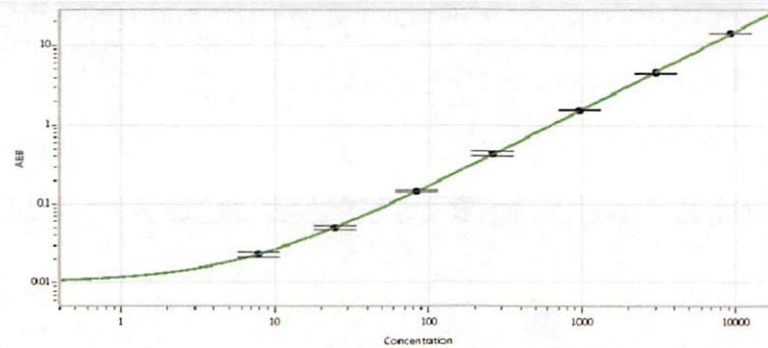
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	9.33
C	29.4
D	104
E	314
F	1165
G	3596
H	9411

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	7.89
C	24.4
D	83.8
E	260
F	971
G	3007
H	9266