

Product Number:	103345
Lot Number:	502892
Expiration:	21-Mar-2022
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103338	111012
Detector Reagent	103339	111207
SBG Reagent	103340	111017
Sample Diluent	103341	111018
Calibrators	103342	112201
Control 1	103343	112202
Control 2	103344	112202
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	112202	2.26	1.73-2.63
Control 2	112202	92.1	73.5-110
Panel 2	111709	3.01	1.47-4.30
Panel 3	111709	28.8	24.6-36.9
Panel 4	111709	83.4	68.2-110

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	112202	8.85	6.73-10.1
Control 2	112202	805	610-915
Panel 2	111709	9.50	7.45-11.2
Panel 3	111709	184	145-218
Panel 4	111709	625	487-731

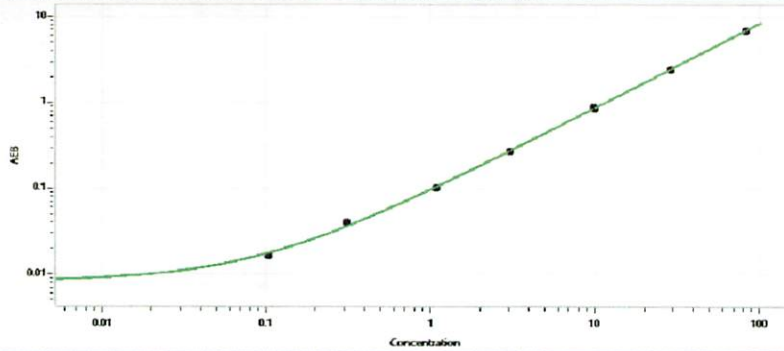
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	112202	150	120-181
Control 2	112202	10849	8199-12298
Panel 2	111709	73.3	62.4-93.8
Panel 3	111709	1453	1186-1778
Panel 4	111709	9614	7766-11952

UCH-L1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	112202	136	101-155
Control 2	112202	11248	8726-13089
Panel 2	111709	112	80.2-135
Panel 3	111709	1790	1417-2126
Panel 4	111709	7621	6242-9363

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

Review/Approval	Junpil Park	QC Supervisor		13 May 2021
	Name	Title	Signature/Date	

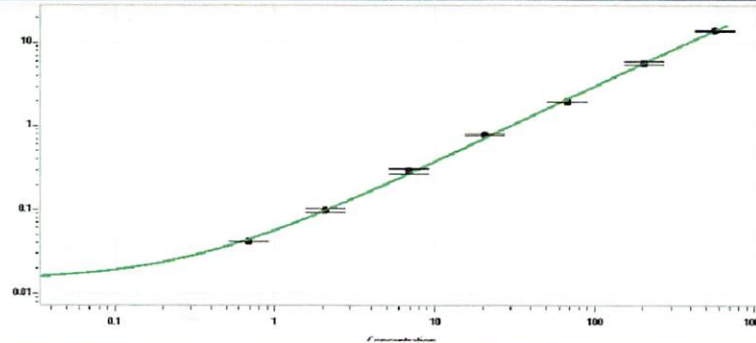
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.096
C	0.297
D	0.957
E	2.84
F	9.12
G	25.8
H	70.9

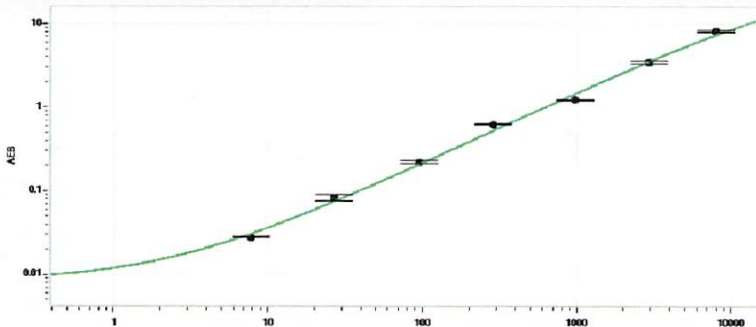
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.646
C	2.05
D	7.02
E	19.8
F	66.9
G	189
H	514

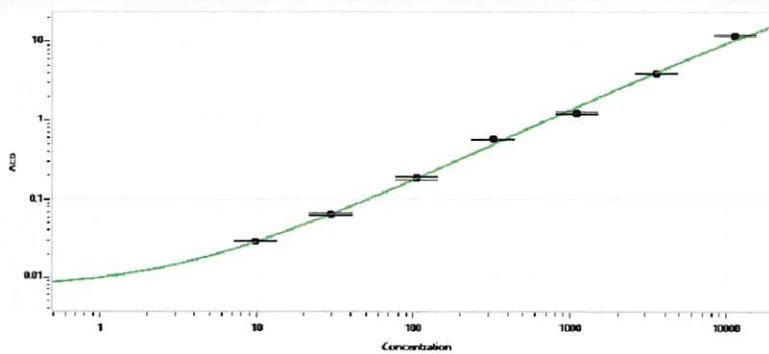
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	9.63
C	30.0
D	103
E	302
F	1056
G	3062
H	8405

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	12.8
C	29.3
D	101
E	305
F	1054
G	3162
H	9438