

Product Number:	102153
Lot Number:	503234
Expiration:	16-Oct-2022
Platform(s):	HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	102141	133203
Detector Reagent	102143	133713
SBG Reagent	102145	132305
Sample Diluent	102147	132306
Calibrators	102150	134311
Control 1	102151	134312
Control 2	102152	134312
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ⁺

¹ SBG Diluent (SR-X) and RGP reagent are not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ Neuro 4-Plex A SR-X Kit Instructions (KI-0031).



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

Release Materials

Tau			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134312	2.69	1.47 - 3.08
Control 2	134312	88.6	71.9 - 108
Panel 1	111709	5.01	1.11 - 7.56
Panel 2	111709	5.32	1.34 - 7.79
Panel 3	111709	25.4	22.5 - 33.7

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134312	4.77	2.69 - 7.55
Control 2	134312	475	370 - 555
Panel 1	111709	4.79	4.12 - 6.23
Panel 2	111709	9.81	7.75 - 11.6
Panel 3	111709	195	147 - 232

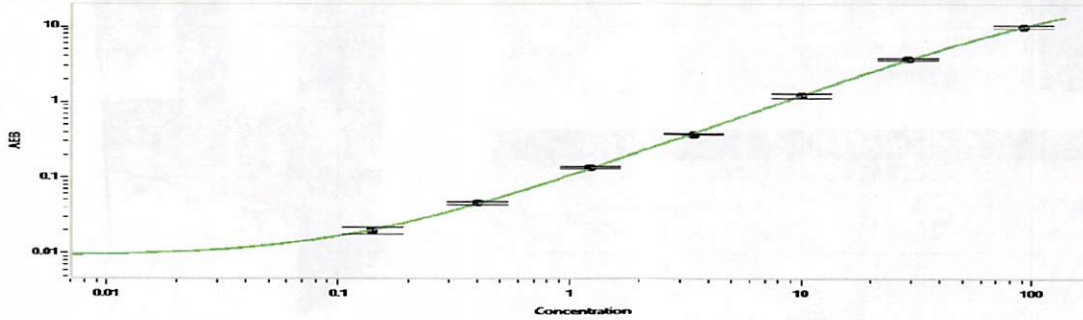
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134312	25.4	17.5 - 39.5
Control 2	134312	1138	878 - 1317
Panel 1	111709	35.3	28.0 - 42.1
Panel 2	111709	108	85.8 - 129
Panel 3	111709	1622	1331 - 1996

UCHL1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134312	111	75.2 - 165
Control 2	134312	9815	7521 - 11281
Panel 1	111709	85.2	69.9 - 123
Panel 2	111709	108	83.9 - 136
Panel 3	111709	1721	1450 - 2176

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

Review/Approval Shivani Goel QC Shivani Goel 06 Jan 2022
Name Title Signature/Date

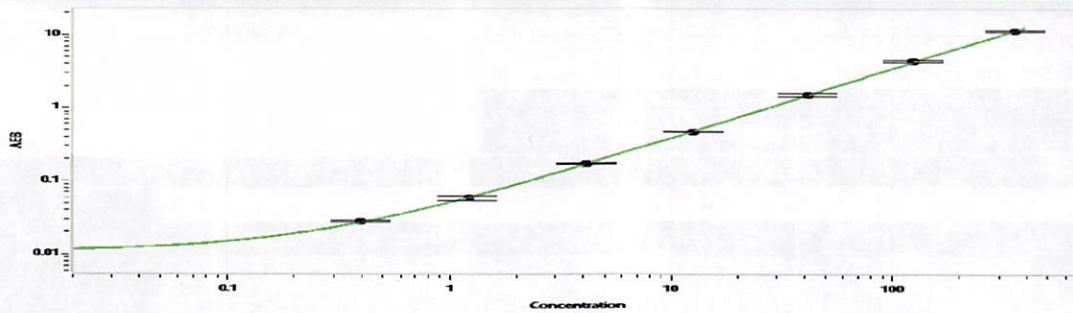
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.140
C	0.400
D	1.23
E	3.44
F	10.0
G	29.0
H	92.1

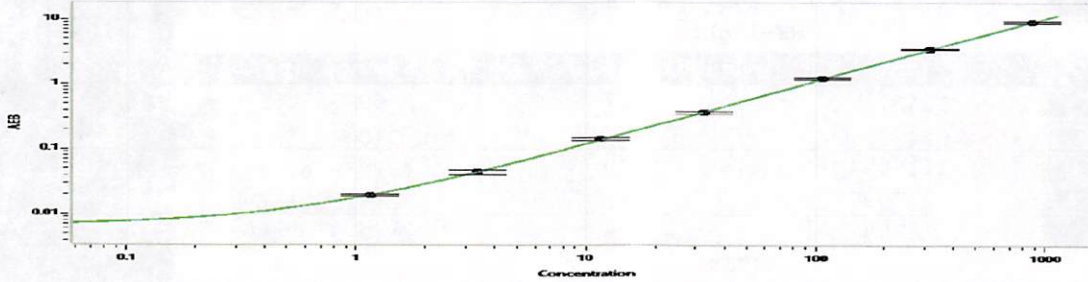
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.396
C	1.21
D	4.15
E	12.6
F	41.2
G	122
H	356

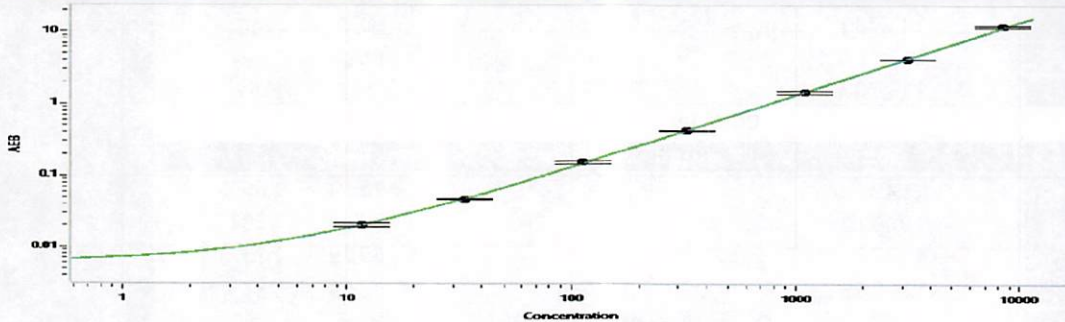
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.15
C	3.36
D	11.5
E	32.8
F	108
G	314
H	872

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	11.6
C	33.3
D	112
E	326
F	1098
G	3136
H	8431