

Product Number:	102153
Lot Number:	502927
Expiration:	22-Apr-2022
Platform(s):	HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	102141	112901
Detector Reagent	102143	113415
SBG Reagent	102145	112907
Sample Diluent	102147	112908
Calibrators	102150	A-E: 113203 F-H: 114804
Control 1	102151	113204
Control 2	102152	113204
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

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).

Release Materials

Tau			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	113204	2.48	1.90 - 2.85
Control 2	113204	99.2	75.2 - 124
Panel 1	111709	2.62	1.73 - 2.94
Panel 2	111709	2.72	2.05 - 3.07
Panel 3	111709	29.9	23.7 - 35.5

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	113204	7.72	6.13 - 9.20
Control 2	113204	744	596 - 894
Panel 1	111709	5.21	4.25 - 6.37
Panel 2	111709	10.0	8.18 - 12.3
Panel 3	111709	190	151 - 226

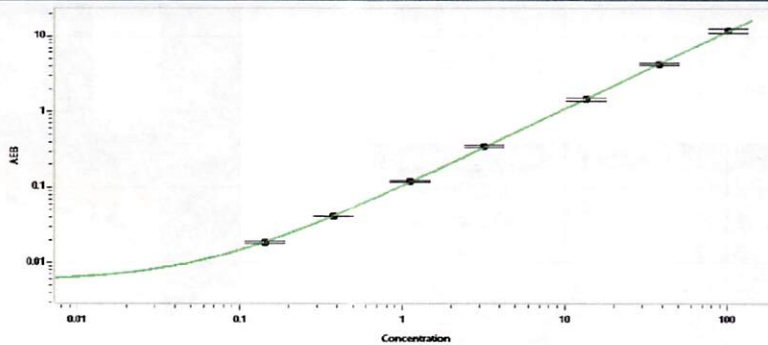
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	113204	32.8	25.1 - 37.6
Control 2	113204	1221	950 - 1524
Panel 1	111709	32.9	25.1 - 37.6
Panel 2	111709	98.8	79.5 - 119
Panel 3	111709	1565	1182 - 1977

UCHL1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	113204	107	81.7 - 140
Control 2	113204	10396	7221 - 14016
Panel 1	111709	88.9	69.0 - 106
Panel 2	111709	110	77.5 - 158
Panel 3	111709	1857	1413-2119

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

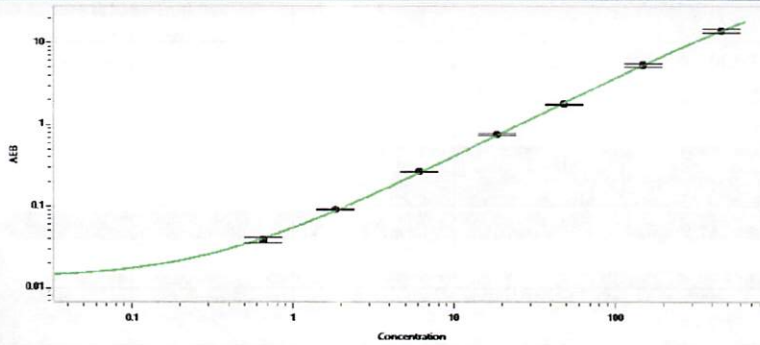
Review/Approval	Junpil Park	QC Supervisor		07 Jun 2021
	Name	Title	Signature/Date	

Tau Calibration Curve



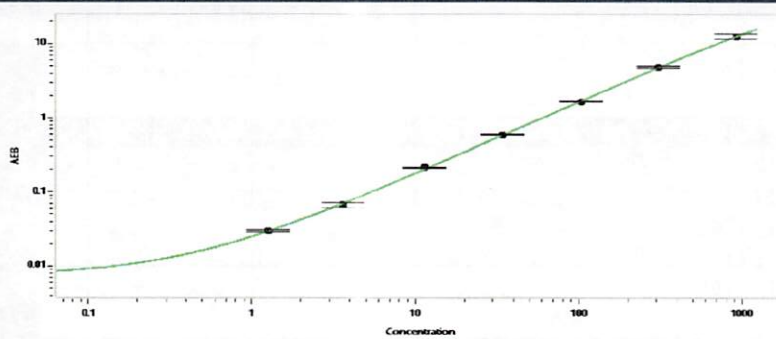
Calibrator Levels (pg/mL)	
A	0.000
B	0.144
C	0.381
D	1.12
E	3.18
F	13.5
G	38.5
H	102

NF-light® Calibration Curve



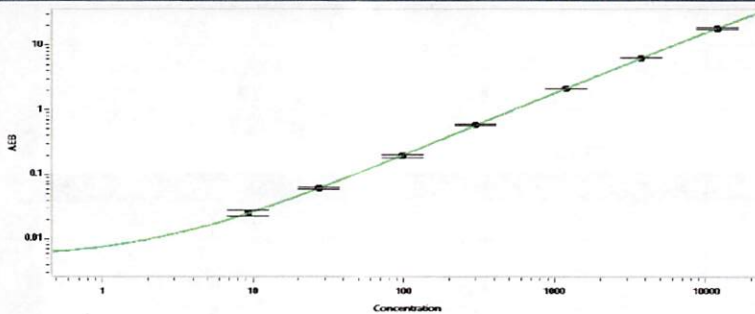
Calibrator Levels (pg/mL)	
A	0.000
B	0.650
C	1.85
D	6.04
E	18.5
F	47.9
G	148
H	455

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.25
C	3.60
D	11.4
E	34.6
F	104
G	305
H	918

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	9.20
C	27.4
D	97.0
E	299
F	1186
G	3696
H	11863