

Product Number:	102153
Lot Number:	503023
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	118009
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	2.70	1.98 - 2.93
Control 2	113204	92.7	75.3 - 121
Panel 1	111709	3.11	1.53 - 3.45
Panel 2	111709	3.47	1.65 - 3.85
Panel 3	111709	30.7	23.8 - 35.8

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	113204	8.37	6.25 - 9.37
Control 2	113204	747	596 - 894
Panel 1	111709	5.64	4.30 - 6.45
Panel 2	111709	11.0	8.31 - 12.5
Panel 3	111709	211	153 - 232

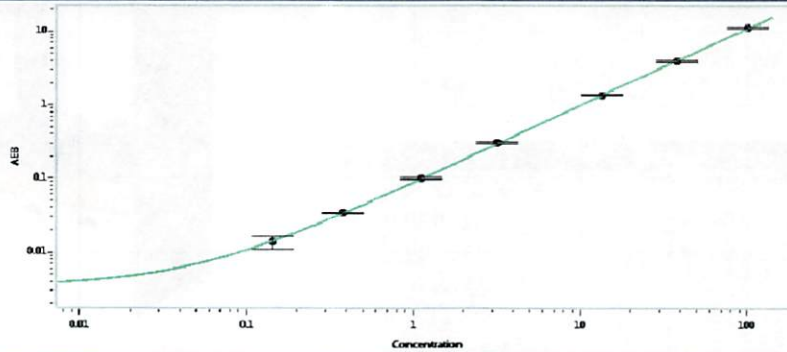
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	113204	33.9	25.5 - 38.2
Control 2	113204	1177	961 - 1489
Panel 1	111709	32.8	25.3 - 38.0
Panel 2	111709	109	81.2 - 122
Panel 3	111709	1716	1212 - 1994

UCHL1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	113204	114	83.2 - 140
Control 2	113204	10060	7407 - 13606
Panel 1	111709	102	64.9 - 116
Panel 2	111709	123	79.1 - 158
Panel 3	111709	2072	1455 - 2182

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

Review/Approval	Junpil Park	QC Supervisor		16 Jul 2021
	Name	Title	Signature/Date	

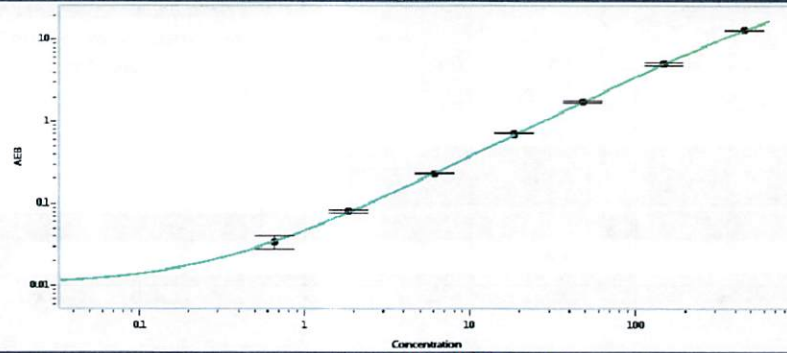
Tau Calibration Curve



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

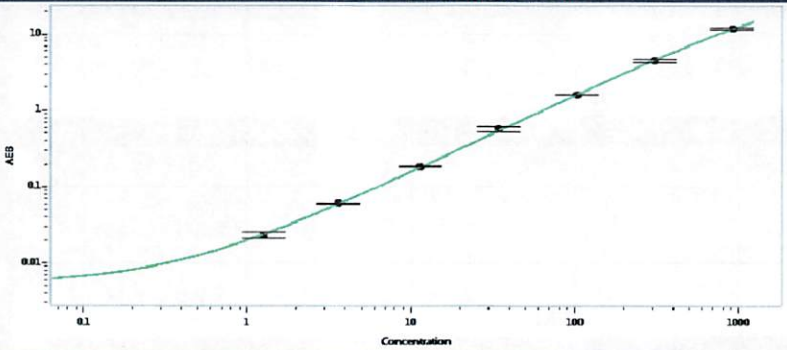
NF-light® Calibration Curve



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

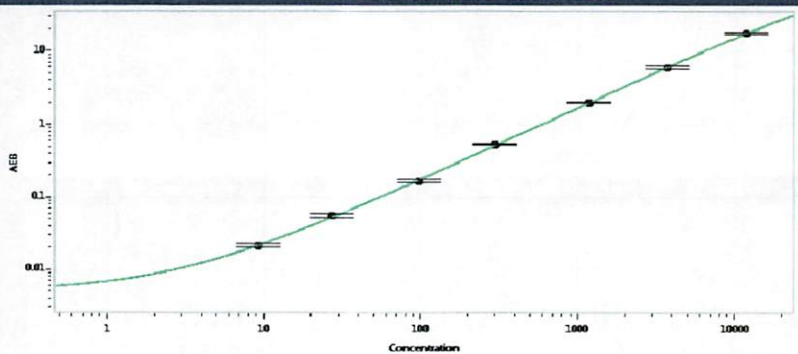
GFAP® Calibration Curve



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