

Product Number:	102153
Lot Number:	502817
Expiration:	9-Dec-2021
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	104813
Detector Reagent	102143	107714
SBG Reagent	102145	034413
Sample Diluent	102147	034418
Calibrators	102150	A-E: 108002 F-H: 108514
Control 1	102151	108003
Control 2	102152	108003
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 A SR-X Kit Instructions (KI-0031).

Release Materials

Tau

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108003	2.36	1.64-2.45
Control 2	108003	90.6	68.2-102
Panel 2	921102	1.67	0.378-1.93
Panel 3	921103	25.7	14.4-38.7

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108003	6.05	4.97-7.50
Control 2	108003	539	443-664
Panel 2	921102	8.27	6.96-10.4
Panel 3	921103	199	151-240


GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108003	29.6	22.4-33.6
Control 2	108003	1142	930-1395
Panel 2	921102	52.1	45.3-68.0
Panel 3	921103	1132	927-1391

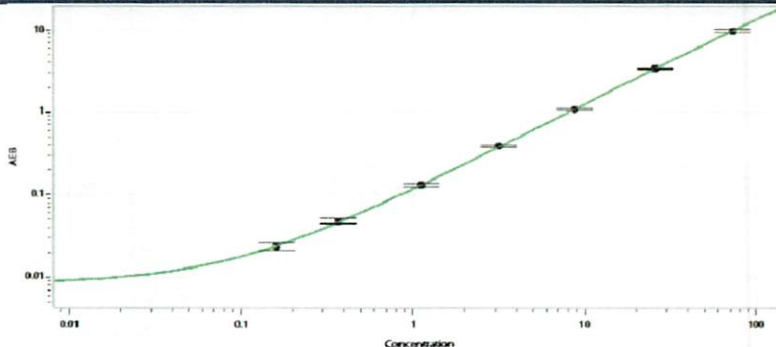
UCHL1®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108003	132	104-155
Control 2	108003	11986	9790-14685
Panel 2	921102	118	57.1-158
Panel 3	921103	3747	1662-5492

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

Review/Approval	Junpil Park	QC Supervisor		31 Mar 2021
	Name	Title	Signature/Date	

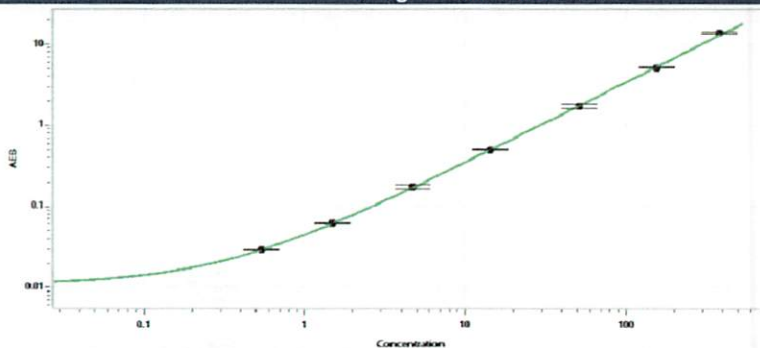
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.162
C	0.370
D	1.12
E	3.17
F	8.76
G	25.7
H	72.8

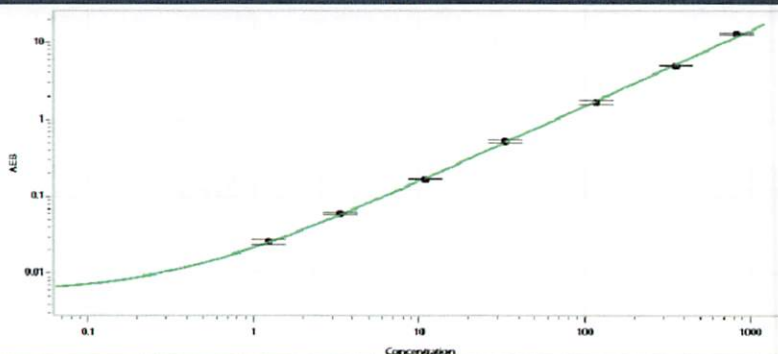
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.542
C	1.49
D	4.70
E	14.3
F	51.5
G	155
H	384

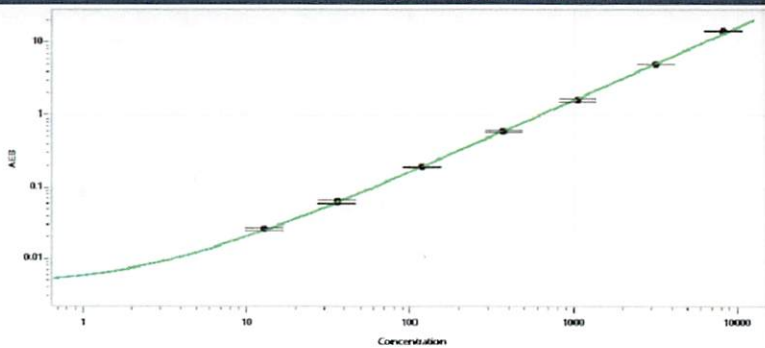
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.24
C	3.34
D	10.9
E	33.4
F	118
G	357
H	829

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	12.9
C	36.0
D	118
E	375
F	1052
G	3168
H	8166