

Product Number:	102153
Lot Number:	503223
Expiration:	16-Oct-2022
Platform(s):	HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	102141	132108
Detector Reagent	102143	132110
SBG Reagent	102145	132109
Sample Diluent	102147	132107
Calibrators	102150	131801
Control 1	102151	131802
Control 2	102152	131802
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹



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

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	131802	2.34	1.71 - 2.85
Control 2	131802	78.5	50.7 - 115
Panel 1	111709	5.00	4.03 - 6.04
Panel 2	111709	4.99	4.03 - 6.24
Panel 3	111709	25.0	22.0 - 33.0

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	131802	5.51	3.98 - 6.72
Control 2	131802	444	214 - 697
Panel 1	111709	4.67	4.02 - 6.03
Panel 2	111709	8.81	7.60 - 11.4
Panel 3	111709	162	131 - 244

GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	131802	23.6	18.4 - 33.3
Control 2	131802	1040	843 - 1473
Panel 1	111709	37.2	28.7 - 43.1
Panel 2	111709	115	89.1 - 134
Panel 3	111709	1566	1357 - 2094

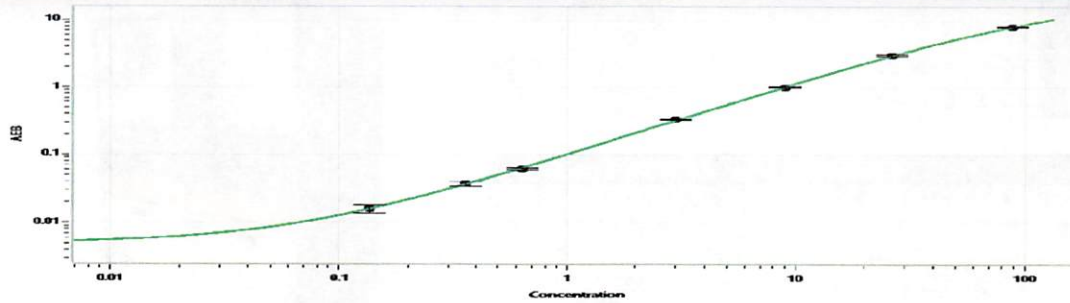
UCHL1®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	131802	133	81.1 - 188
Control 2	131802	9374	7781 - 11856
Panel 1	111709	98.3	76.5 - 127
Panel 2	111709	107	76.8 - 142
Panel 3	111709	1663	1455 - 2182

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

Review/Approval: Junpil Park (Name), QC Supervisor (Title), [Signature] 29 NOV 2021 (Signature/Date)

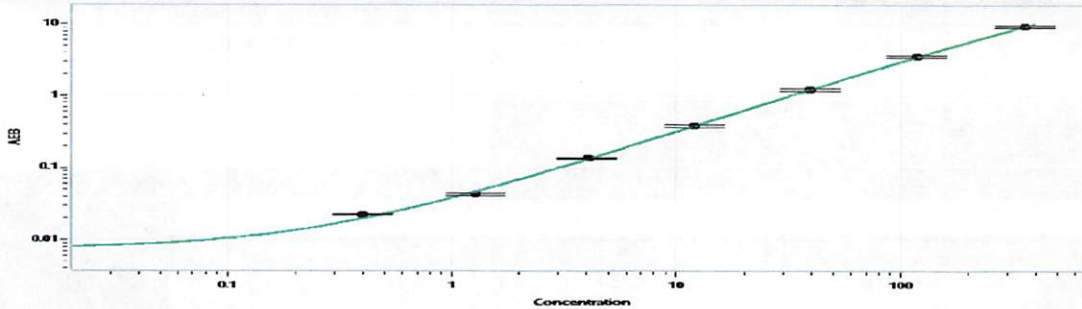
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.136
C	0.360
D	0.636
E	2.98
F	9.03
G	26.3
H	88.7

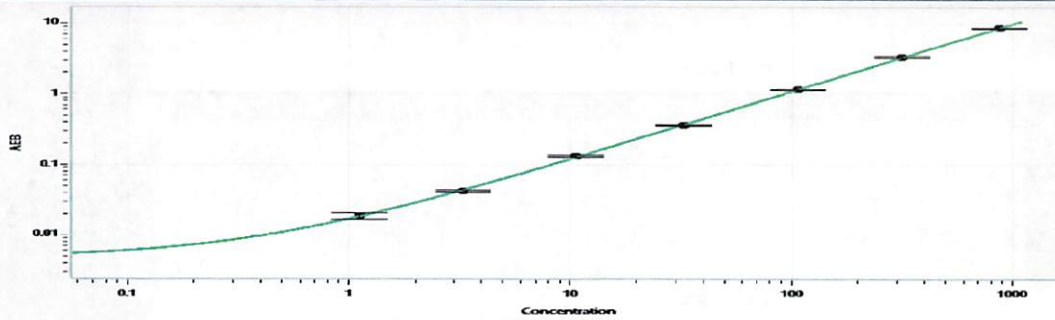
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.400
C	1.28
D	4.01
E	12.1
F	39.6
G	117
H	359

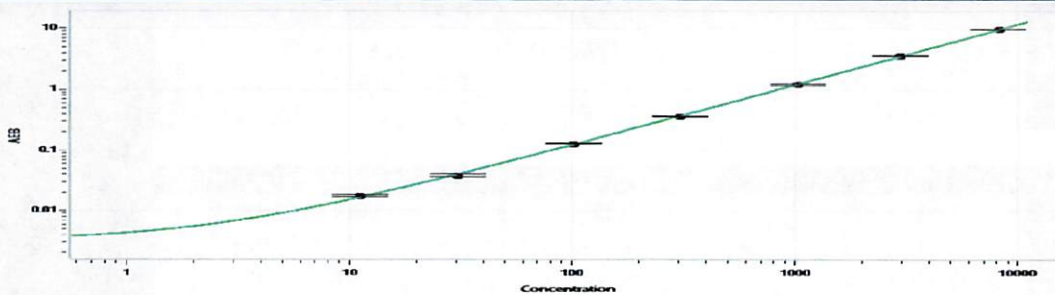
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.11
C	3.29
D	10.6
E	32.5
F	107
G	316
H	866

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	11.1
C	30.5
D	100
E	305
F	1021
G	2932
H	8156