

Product Number:	102153
Lot Number:	502509
Expiration:	3-Jun-2021
Platform(s):	HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	102141	023404
Detector Reagent	102143	020509
SBG Reagent	102145	017802
Sample Diluent	102147	017805
Calibrators	102150	019901
Control 1	102151	019902
Control 2	102152	019902
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	019902	2.17	1.66-2.50
Control 2	019902	87.2	70.1-105
Panel 1	921102	0.914	0.311-1.73
Panel 2	921103	0.757	14.4-38.7
Panel 3	921104	24.3	27.7-53.0

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	019902	6.29	5.13-7.70
Control 2	019902	504	428-642
Panel 1	921102	9.90	7.00-10.5
Panel 2	921103	7.36	151-240
Panel 3	921104	177	561-897

GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	019902	21.6	16.7-27.4
Control 2	019902	992	832-1248
Panel 1	921102	76.3	45.7-68.5
Panel 2	921103	52.2	927-1391
Panel 3	921104	969	3970-8978

UCHL1®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	019902	90.5	82.0-127
Control 2	019902	8049	7183-10774
Panel 1	921102	119	52.9-161
Panel 2	921103	116	1662-5492
Panel 3	921104	3471	4277-15688

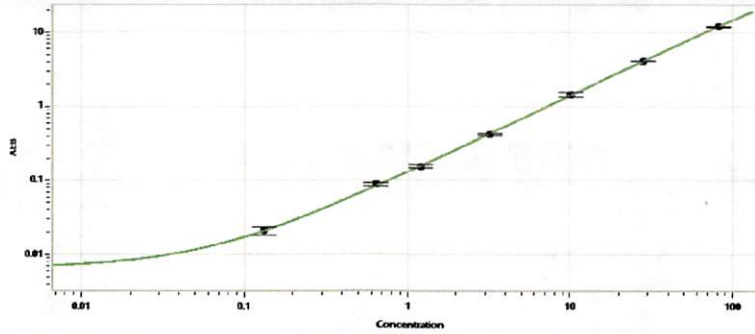
Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.



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

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 23Sep2020
	Name	Title	Signature/Date

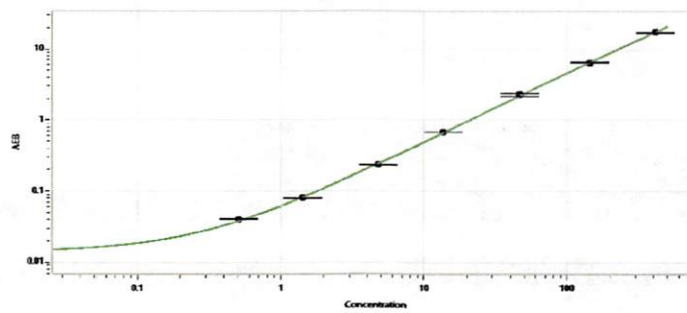
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.133
C	0.647
D	1.22
E	3.22
F	10.1
G	27.9
H	81.1

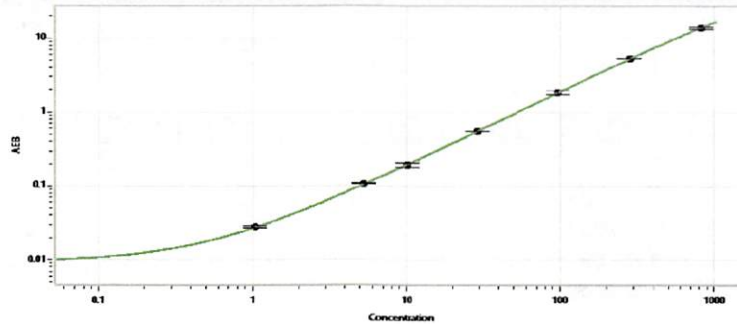
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.506
C	1.43
D	4.81
E	13.7
F	46.4
G	144
H	410

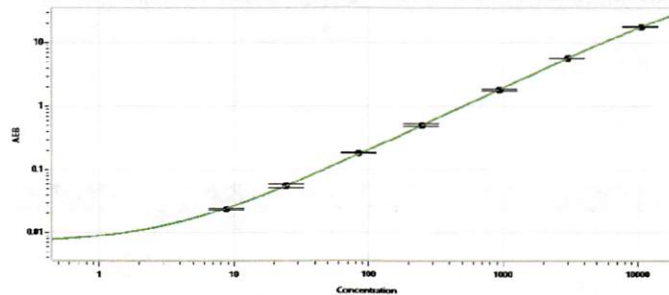
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.04
C	5.29
D	10.1
E	28.8
F	97.0
G	282
H	820

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	8.77
C	24.4
D	84.4
E	248
F	940
G	2998
H	10508