

Product Number:	102153
Lot Number:	502851
Expiration:	7-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	109705
Detector Reagent	102143	111819
SBG Reagent	102145	109813
Sample Diluent	102147	109814
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	2.41	1.60-2.71
Control 2	108003	92.7	70.0-105
Panel 2	921102	1.36	0.683-1.82
Panel 3	921103	25.3	14.4-38.7
Panel 4	921104	39.5	27.7-53.0

NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108003	6.49	5.05-7.58
Control 2	108003	551	442-663
Panel 2	921102	9.60	6.96-10.4
Panel 3	921103	188	151-240
Panel 4	921104	667	561-897

GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108003	31.4	23.0-35.1
Control 2	108003	1171	932-1398
Panel 2	921102	55.3	45.3-68.0
Panel 3	921103	975	927-1391
Panel 4	921104	4536	3970-8978

UCHL1®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108003	136	105-158
Control 2	108003	12160	9771-14657
Panel 2	921102	120	57.1-158
Panel 3	921103	3472	1662-5492
Panel 4	921104	8062	4277-15688

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

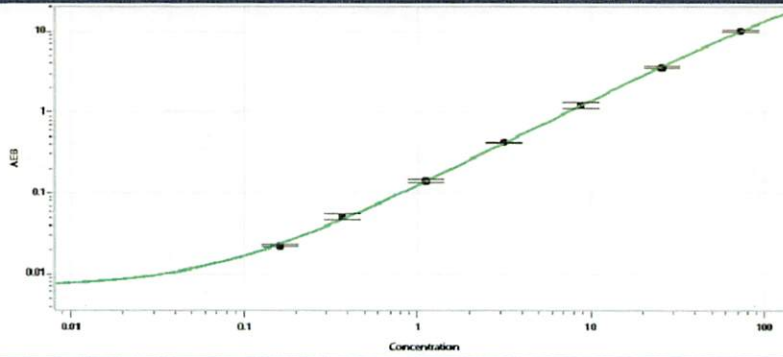
Review/Approval

Junpil Park
Name

QC Supervisor
Title

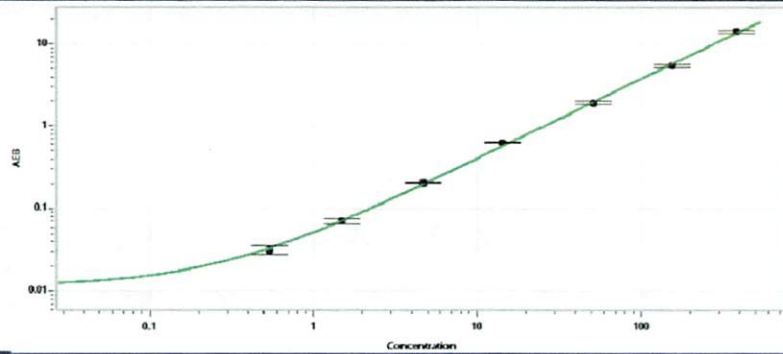
 04 May 2021
Signature/Date

Tau Calibration Curve



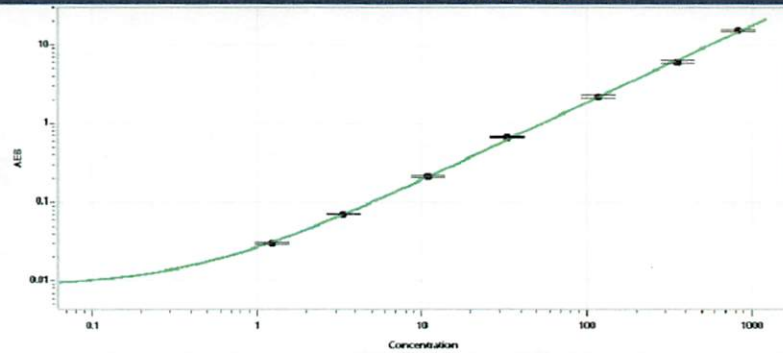
Calibrator Levels (pg/mL)	
A	0.000
B	0.162
C	0.370
D	1.12
E	3.17
F	8.76
G	25.7
H	72.8

NF-light® Calibration Curve



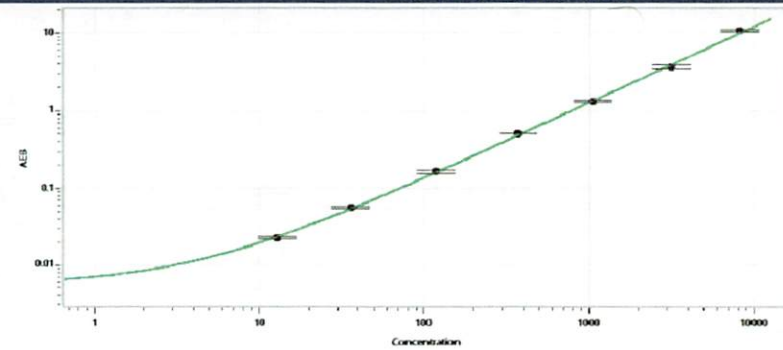
Calibrator Levels (pg/mL)	
A	0.000
B	0.542
C	1.49
D	4.70
E	14.3
F	51.5
G	155
H	384

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