

Product Number:	102153
Lot Number:	503478
Expiration:	4-May-2023
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	216006
Detector Reagent	102143	225710
SBG Reagent	102145	216103
Sample Diluent	102147	216205
Calibrators	102150	215703
Control 1	102151	215704
Control 2	102152	215704
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)	Spec. Range (pg/mL)
Control 1	215704	2.15	1.74 - 2.61
Control 2	215704	79.6	62.9 - 106
Panel 1	206807	5.92	4.36 - 13.3
Panel 2	206807	5.18	4.08 - 7.93
Panel 3	206807	20.8	16.5 - 25.9

NF-light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	215704	4.06	3.01 - 5.11
Control 2	215704	405	333 - 500
Panel 1	206807	10.7	6.25 - 12.5
Panel 2	206807	4.32	1.77 - 5.31
Panel 3	206807	147	91.9 - 161

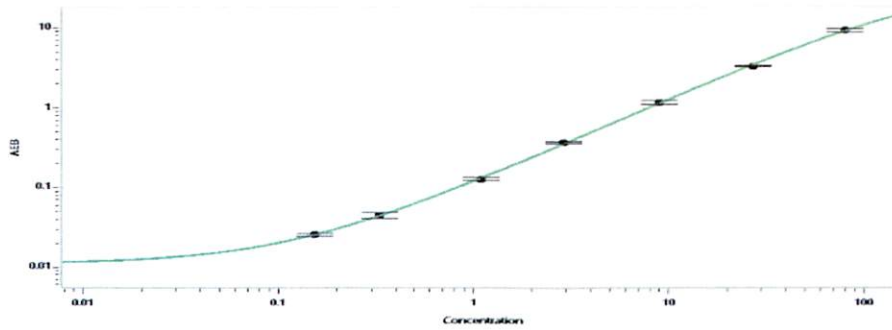
GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	215704	21.9	17.2 - 25.7
Control 2	215704	1109	893 - 1340
Panel 1	206807	62.6	55.6 - 83.4
Panel 2	206807	23.9	21.1 - 31.7
Panel 3	206807	816	680 - 1026

UCHL1®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	215704	141	109 - 172
Control 2	215704	14075	11094 - 16641
Panel 1	206807	94.7	43.8 - 117
Panel 2	206807	85.9	48.9 - 109
Panel 3	206807	1234	916 - 1405

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

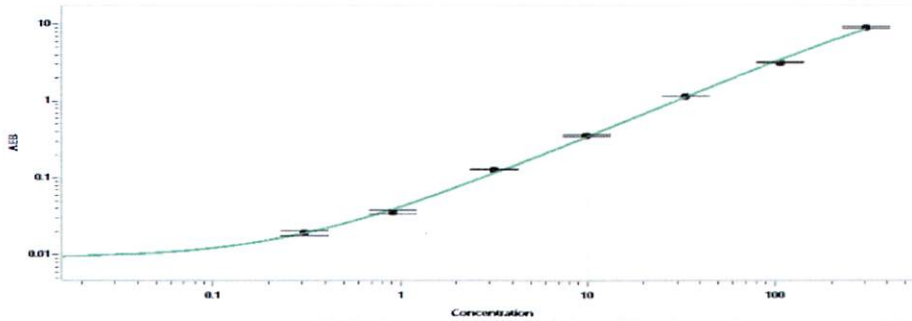
Review/Approval	<u>Sonia Khillan</u> Name	QC Title	<u>Sonia Khillan / 27SEP2022</u> Signature/Date
-----------------	------------------------------	-------------	--

Tau Calibration Curve



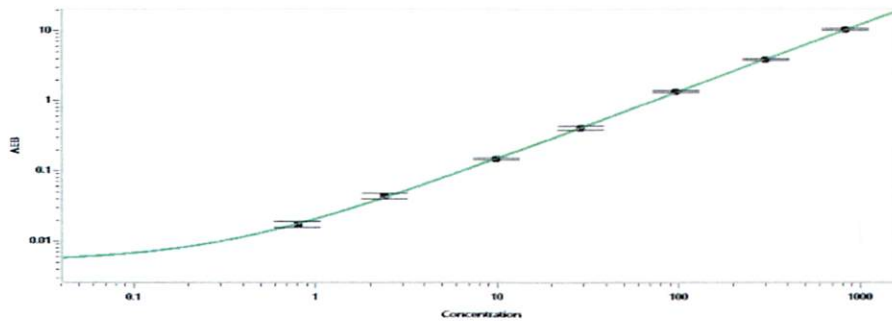
Calibrator Levels (pg/mL)	
A	0
B	0.155
C	0.331
D	1.09
E	2.93
F	8.95
G	26.9
H	79.7

NF-light® Calibration Curve



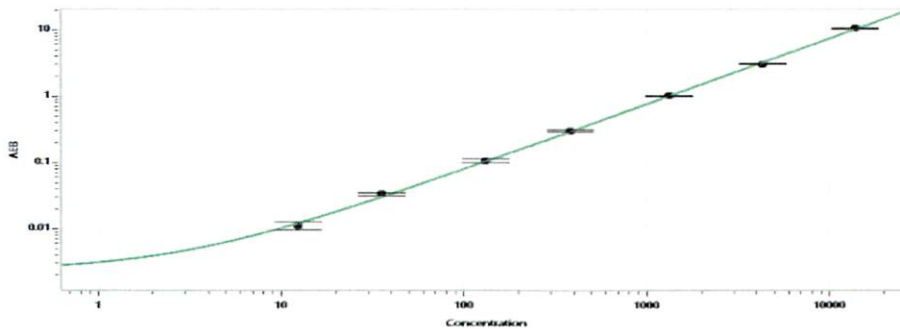
Calibrator Levels (pg/mL)	
A	0
B	0.311
C	0.918
D	3.16
E	9.85
F	33.2
G	107
H	305

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0
B	0.848
C	2.43
D	9.88
E	28.7
F	96.2
G	301
H	827

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)	
A	0
B	12.5
C	35.6
D	131
E	382
F	1335
G	4329
H	13889