

Product Number:	102153
Lot Number:	503291
Expiration:	13-Jan-2023
Platform(s):	HD-X & SR-X



Data below represents results generated on the Simoa® HD-X / SR-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102141	201305
Detector Reagent	102143	201306
SBG Reagent	102145	201307
Sample Diluent	102147	201308
Calibrators	102150	209406
Control 1	102151	209407
Control 2	102152	209407
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)	Mean Range (pg/mL)
Control 1	209407	2.10	1.65-2.48
Control 2	209407	64.6	55.1-82.7
Panel 1	206807	11.2	6.77-12.4
Panel 2	206807	7.00	4.44-8.02
Panel 3	206807	22.9	16.1-26.4

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	209407	4.81	3.36-6.05
Control 2	209407	380	326-489
Panel 1	206807	9.81	6.18-11.9
Panel 2	206807	3.98	1.71-4.94
Panel 3	206807	126	97.2-146

GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	209407	20.9	15.4-27.3
Control 2	209407	944	790-1186
Panel 1	206807	71.5	56.9-85.4
Panel 2	206807	27.2	21.7-32.6
Panel 3	206807	959	690-1035

UCHL1®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	209407	136	103-166
Control 2	209407	11713	9969-14954
Panel 1	206807	88.9	49.2-105
Panel 2	206807	79.5	46.3-109
Panel 3	206807	1272	902-1383

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

Review/Approval

Shivani Goel

QC Manager

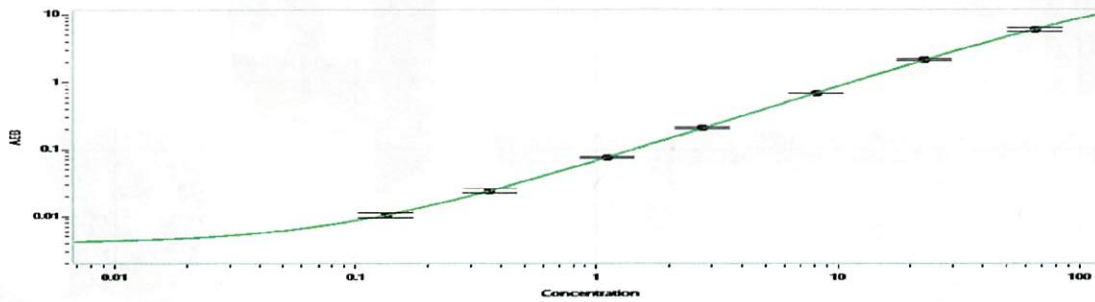
Shivani Goel 12 May 2022

Name

Title

Signature/Date

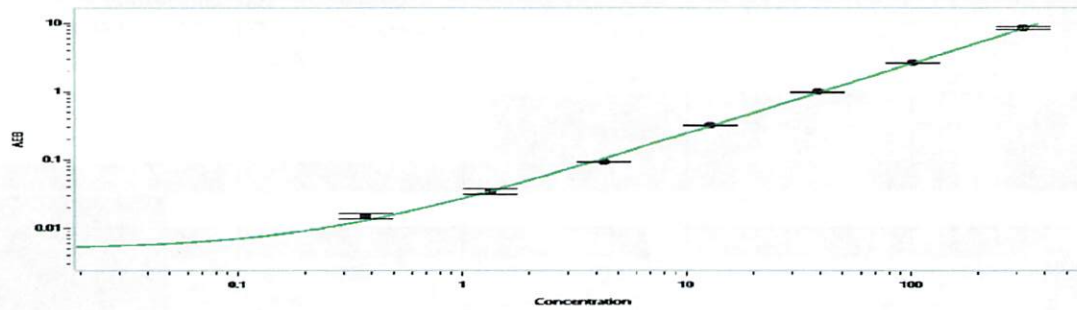
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.134
C	0.360
D	1.11
E	2.75
F	8.15
G	22.7
H	65.4

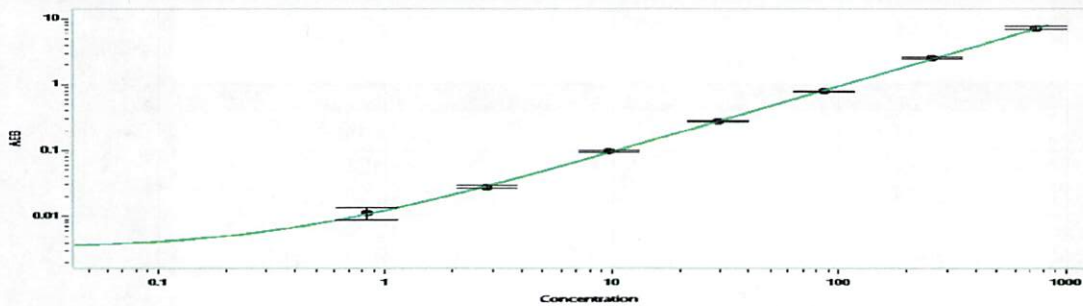
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.373
C	1.35
D	4.34
E	12.8
F	38.7
G	102
H	319

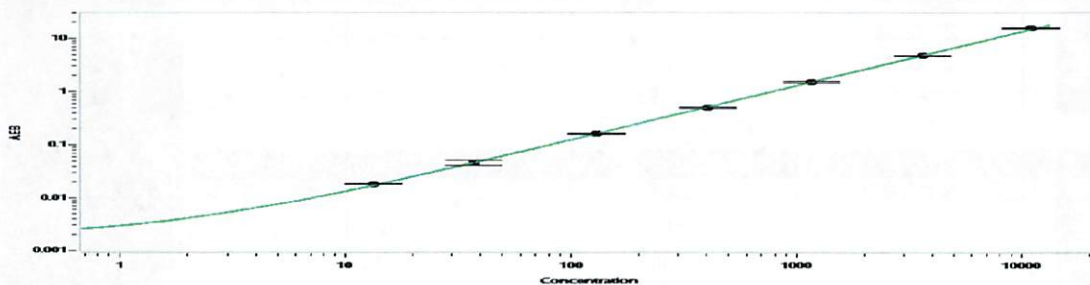
GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.841
C	2.84
D	9.78
E	29.6
F	86.7
G	257
H	744

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	13.3
C	37.5
D	131
E	404
F	1164
G	3616
H	10975