

Product Number:	102153
Lot Number:	503770
Expiration:	3-Mar-2024
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	306204
Detector Reagent	102143	306514
SBG Reagent	102145	307410
Sample Diluent	102147	306515
Calibrators	102150	314001 ²
Control 1	102151	314002
Control 2	102152	314002
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

² Calibrators Control Lot number is 314001A

† 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	314002	2.02	1.63 - 2.45
Control 2	314002	68.2	53.8 - 80.7
Panel 1	308106	28.1	25.7 - 39.0
Panel 2	309707	2.65	2.05 - 3.17
Panel 3	311701	73.4	60.1 - 90.2

NF-light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	314002	5.78	4.53 - 6.80
Control 2	314002	610	447 - 670
Panel 1	308106	11.1	8.85 - 13.3
Panel 2	309707	7.77	6.28 - 9.42
Panel 3	311701	496	406 - 608

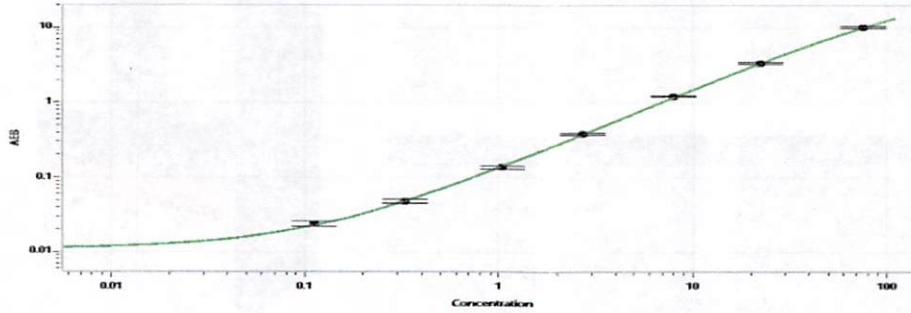
GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	314002	20.3	16.0 - 24.0
Control 2	314002	1057	783 - 1174
Panel 1	308106	48.7	31.7 - 52.9
Panel 2	309707	85.4	64.2 - 96.3
Panel 3	311701	1151	883 - 1324

UCHL1®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	314002	132	108 - 163
Control 2	314002	11301	9084 - 13626
Panel 1	308106	88.9	54.2 - 108
Panel 2	309707	100	83.4 - 132
Panel 3	311701	6319	5551 - 8326

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

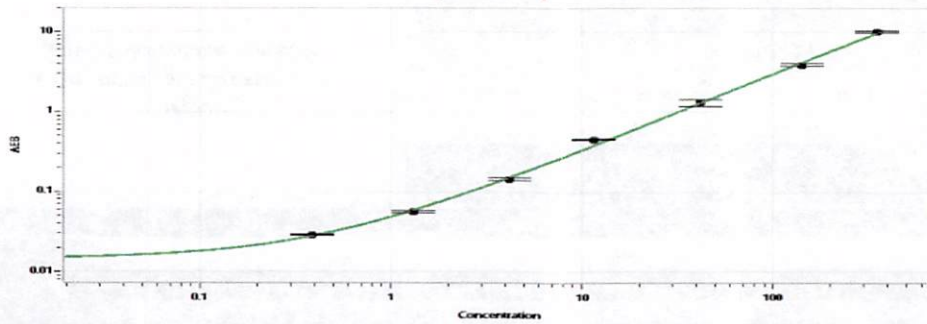
Review/Approval Sonia Khillan QC Sonia Khillan 06 JUN 2023
Name Title Signature/Date

Tau Calibration Curve



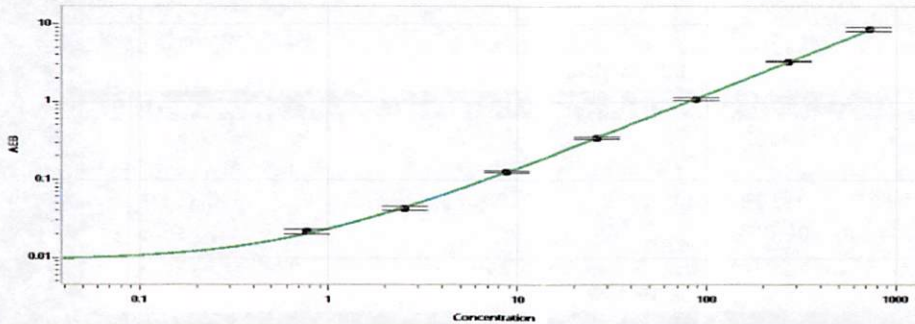
Calibrator Levels (pg/mL)	
A	0.000
B	0.112
C	0.330
D	1.04
E	2.70
F	8.00
G	22.4
H	75.7

NF-light® Calibration Curve



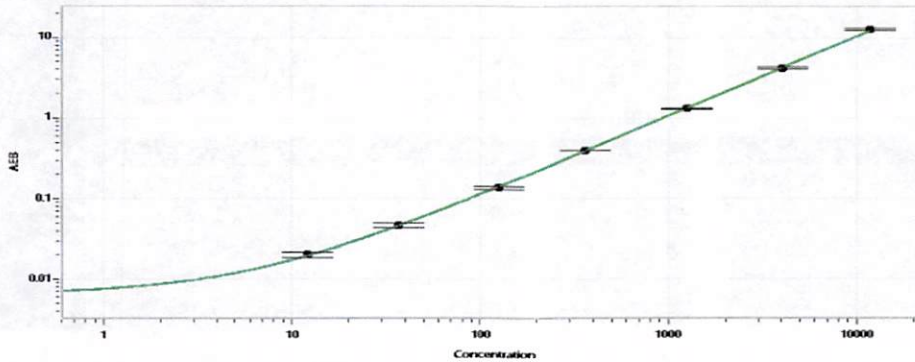
Calibrator Levels (pg/mL)	
A	0.000
B	0.386
C	1.31
D	4.19
E	11.5
F	41.9
G	142
H	352

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.771
C	2.55
D	8.71
E	26.3
F	88.3
G	273
H	723

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	12.0
C	37.0
D	126
E	362
F	1252
G	4022
H	11702