

Product Number:	102153
Lot Number:	503676
Expiration:	30-Nov-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	233404
Detector Reagent	102143	233505
SBG Reagent	102145	233506
Sample Diluent	102147	233507
Calibrators	102150	232013
Control 1	102151	232014
Control 2	102152	233014
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	232014	2.30	1.93 - 2.89
Control 2	232014	83.9	69.2 - 121
Panel 1	206807	5.14	2.90 - 13.6
Panel 2	206807	3.80	2.88 - 8.36
Panel 3	206807	18.3	15.8 - 25.7

NF-light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	232014	4.85	3.96 - 7.19
Control 2	232014	510	405 - 657
Panel 1	206807	8.36	6.16 - 12.3
Panel 2	206807	3.14	1.81 - 5.13
Panel 3	206807	149	91.2 - 169

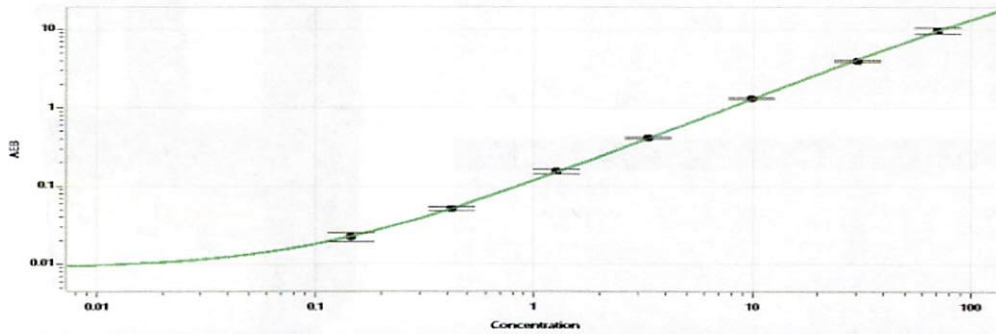
GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	232014	20.5	17.5 - 26.3
Control 2	232014	1122	901 - 1471
Panel 1	206807	57.4	51.9 - 83.0
Panel 2	206807	20.3	18.1 - 32.6
Panel 3	206807	958	676 - 1065

UCHL1®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	232014	135	100 - 156
Control 2	232014	11048	8149 - 14778
Panel 1	206807	94.3	46.5 - 120
Panel 2	206807	89.6	50.9 - 110
Panel 3	206807	1432	864 - 1546

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

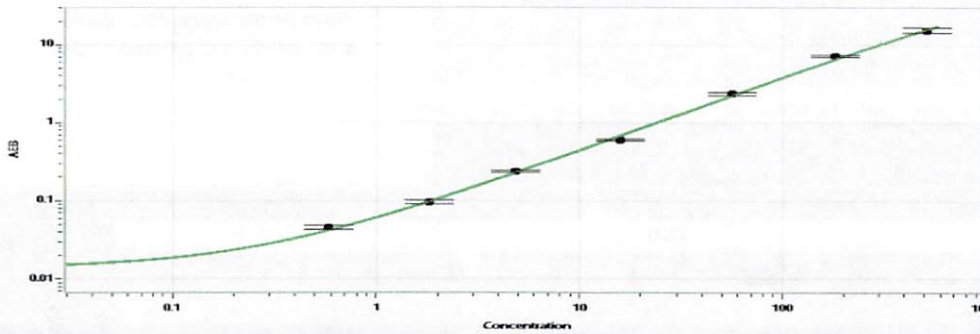
Review/Approval Sonia Khillan QC Sonia Khillan 09NOV2023
Name Title Signature/Date

Tau Calibration Curve



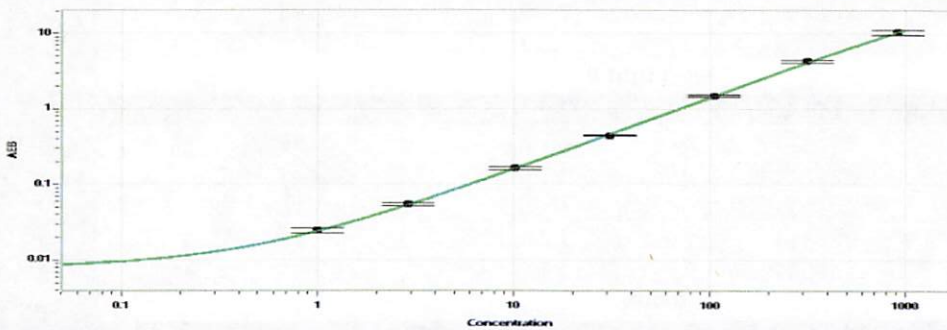
Calibrator Levels (pg/mL)	
A	0.00
B	0.145
C	0.421
D	1.26
E	3.32
F	9.95
G	30.1
H	93.8

NF-light® Calibration Curve



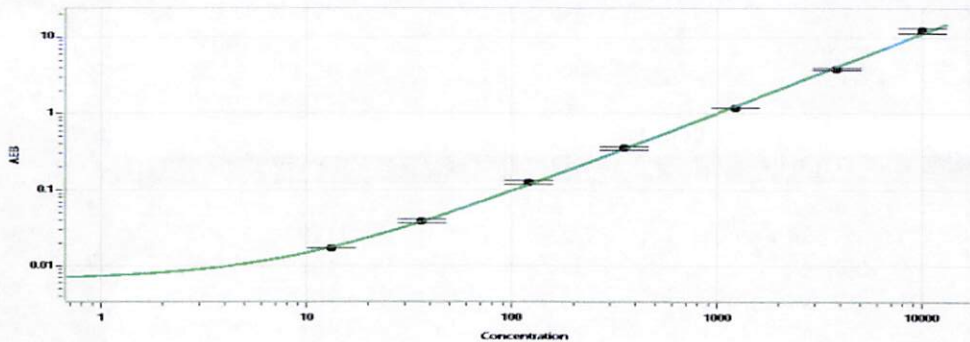
Calibrator Levels (pg/mL)	
A	0.00
B	0.589
C	1.83
D	4.84
E	15.9
F	55.9
G	181
H	515

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.00
B	1.00
C	2.93
D	10.2
E	31.0
F	106
G	316
H	920

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.00
B	13.3
C	36.0
D	120
E	349
F	1220
G	3784
H	13149