

Product Number:	102153
Lot Number:	503857
Expiration:	19-May-2024
Platform(s):	HD-X & SR-X



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

Component	Part Number	Lot Number	Control Number
Bead Reagent	102141	313904	313904A
Detector Reagent	102143	317711	317711A
SBG Reagent	102145	314305	314305A
Sample Diluent	102147	314306	314306A
Calibrators	102150	320603	320603
Control 1	102151	320604	320604
Control 2	102152	320604	320604
SBG Diluent (SR-X)†	103158	N/A ¹	N/A ¹
RGP Reagent	103159	N/A ¹	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	320604	2.46	1.95 - 2.93
Control 2	320604	79.6	60.3 - 108
Panel 1	308106	30.0	25.6 - 38.4
Panel 2	309707	2.91	2.05 - 3.29
Panel 3	311701	71.7	59.5 - 89.3

NF-light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	320604	5.60	4.43 - 6.93
Control 2	320604	497	382 - 705
Panel 1	308106	9.00	8.03 - 13.3
Panel 2	309707	6.87	6.06 - 9.25
Panel 3	311701	417	388 - 590

GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	320604	26.1	19.1 - 28.6
Control 2	320604	1183	912 - 1368
Panel 1	308106	45.7	32.7 - 53.2
Panel 2	309707	79.1	64.0 - 96.0
Panel 3	311701	1113	884 - 1326

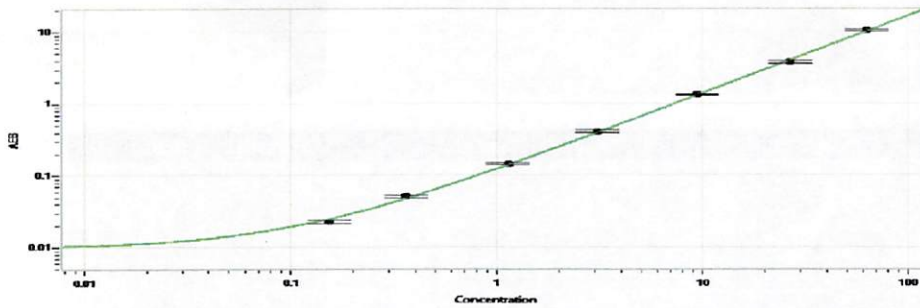
UCHL1®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	320604	150	114 - 178
Control 2	320604	12287	8430 - 17010
Panel 1	308106	122	31.2 - 150
Panel 2	309707	120	81.7 - 142
Panel 3	311701	6385	5407 - 8145

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

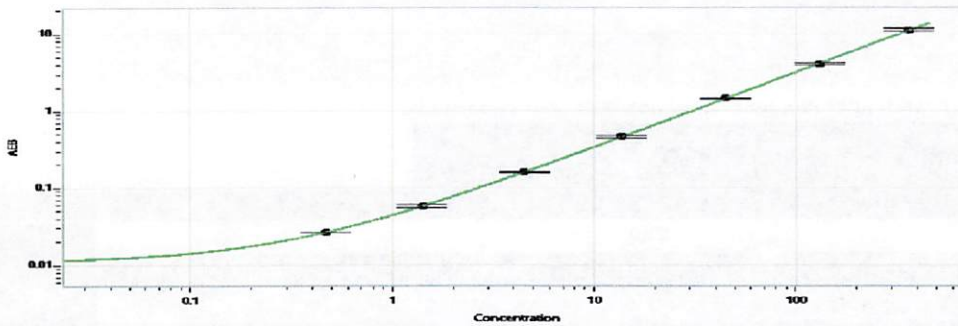
Review/Approval Sonia Khilka QC Sonia Khilka 09NOV2023
 Name Title Signature/Date

Tau Calibration Curve



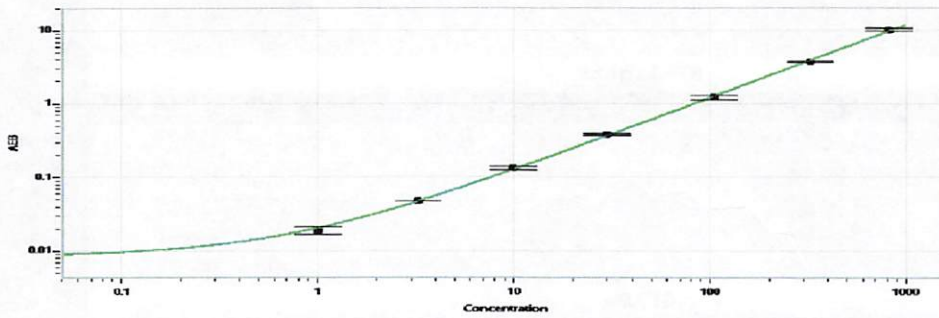
Calibrator Levels (pg/mL)	
A	0.000
B	0.154
C	0.363
D	1.13
E	3.06
F	9.44
G	26.4
H	77.9

NF-light® Calibration Curve



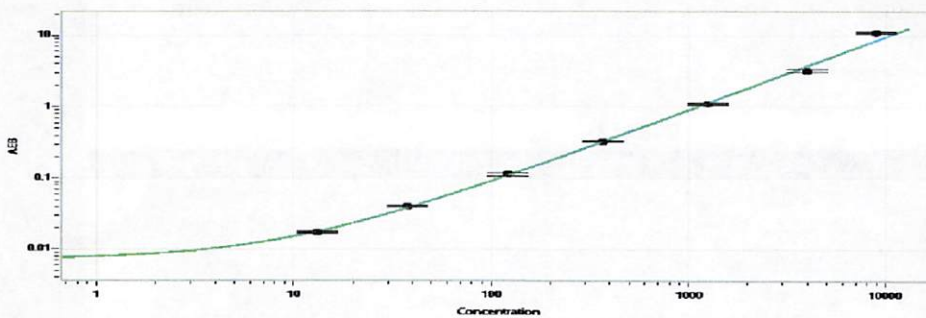
Calibrator Levels (pg/mL)	
A	0.000
B	0.466
C	1.41
D	4.52
E	13.6
F	45.1
G	132
H	362

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.00
C	3.28
D	10.0
E	29.9
F	105
G	322
H	815

UCHL1® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	13.2
C	37.3
D	121
E	366
F	1248
G	3963
H	11889