

Product Number:	104080
Lot Number:	503878
Expiration:	26-Apr-2024
Platform(s):	HD-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	103510	311608	311608A
Detector Reagent	103512	311802	311802A
SBG Reagent	103514	311803	311803A
Sample Diluent	103516	311804	311804A
Calibrators	103517	313004	313004A
Control 0	104079	313104	313104A
Control 1	104078	313104	313104A
Control 2	104077	313104	313104A
RGP Reagent	103159	N/A ¹	N/A ¹

¹ RGP / SBG Diluent (SR-X) are not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa® Neuro 2-Plex B SR-X Kit Instructions (KI-0062).

Release Materials

NF-light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 0	313104	7.99	6.76 - 10.14
Control 1	313104	136	102 - 152
Control 2	313104	407	307 - 460
Panel 1	310007	8.21	4.41 - 9.34
Panel 2	310007	5.31	2.27 - 6.19
Panel 3	306904	107	73.1 - 110
Panel 4	306904	527	353 - 584

GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 0	313104	71.1	50.3 - 81.1
Control 1	313104	411	321 - 555
Control 2	313104	1988	1714 - 2572
Panel 1	310007	68.6	43.7 - 86.7
Panel 2	310007	63.9	39.9 - 71.8
Panel 3	306904	1177	775 - 1475
Panel 4	306904	7849	6304 - 9456

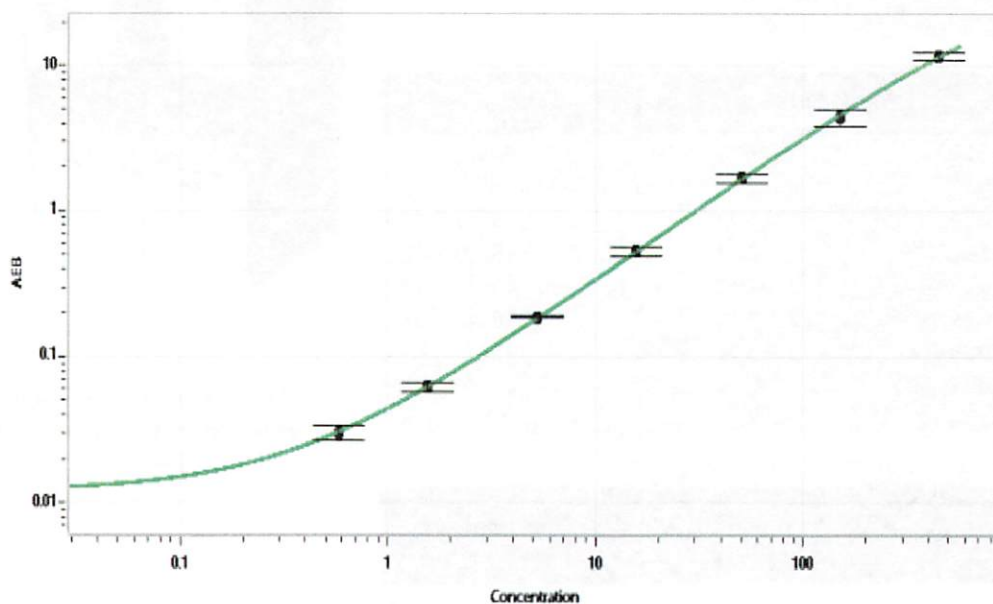
Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Review/Approval

Sonia Khillan QC
Name Title

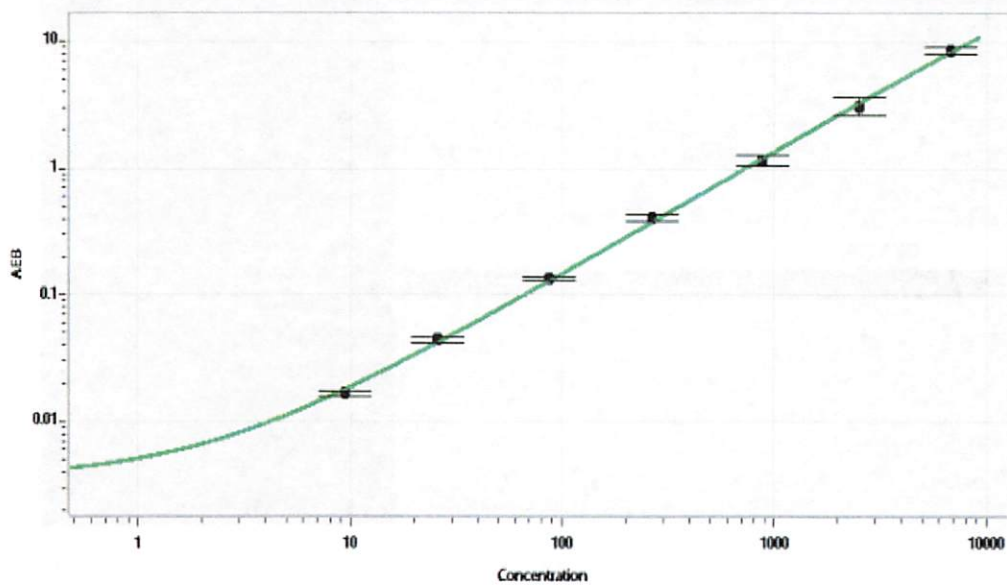
[Signature] 06 JUN 2023
Signature/Date

NF-light® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.579
C	1.56
D	5.21
E	15.5
F	50.5
G	149
H	444

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	9.41
C	25.8
D	86.5
E	267
F	879
G	2507
H	6823