

Product Number:	104080
Lot Number:	503921
Expiration:	21-Jun-2024
Platform(s):	HD-X



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

Component	Part Number	Lot Number	Control Number
Bead Reagent	103510	317211	317211A
Detector Reagent	103512	321313	N/A
SBG Reagent	103514	320905	N/A
Sample Diluent	103516	317307	317307A
Calibrators	103517	321906	N/A
Control 0	104079	319105	319105A
Control 1	104078	319105	319105A
Control 2	104077	319105	319105A
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	319105	3.30	2.54 - 5.87
Control 1	319105	72.7	61.3 - 92.0
Control 2	319105	219	185 - 277
Panel 1	310007	6.53	5.26 - 11.1
Panel 2	323304	5.56	4.46 - 8.31
Panel 3	306904	85.3	74.0 - 111
Panel 4	306904	466	359 - 597

GFAP®

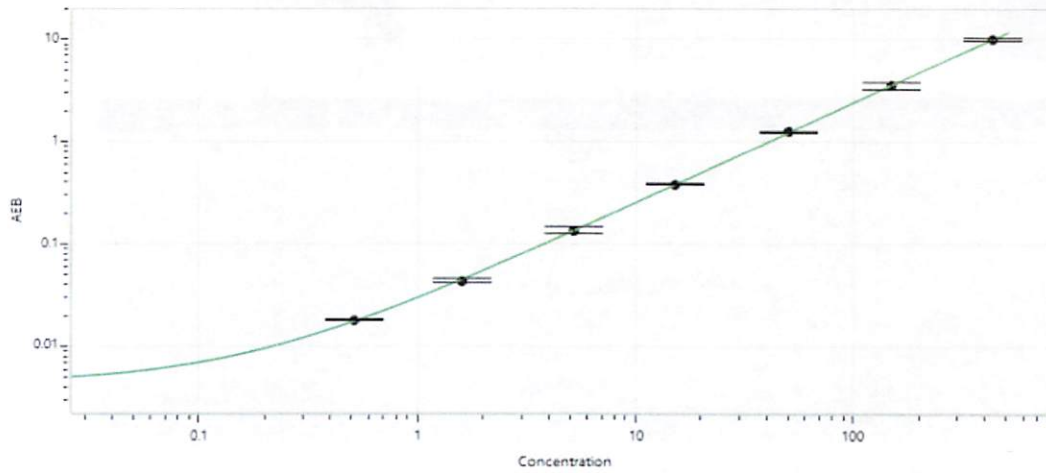
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 0	319105	73.7	41.6 - 79.0
Control 1	319105	456	309 - 464
Control 2	319105	2258	1561 - 2341
Panel 1	310007	85.7	63.0 - 105
Panel 2	323304	66.3	34.3 - 82.1
Panel 3	306904	1314	795 - 1440
Panel 4	306904	8508	6199 - 9299

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

Review/Approval

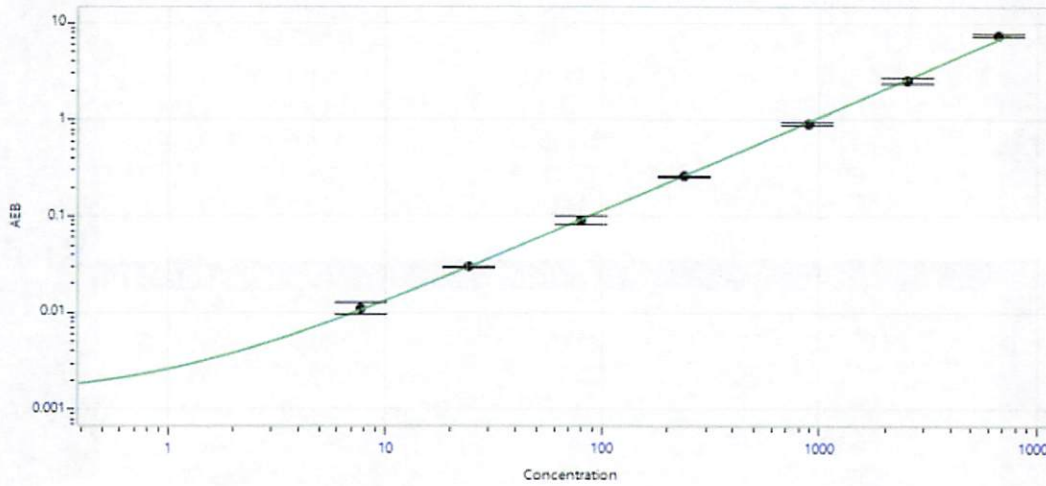
Sonia Khillan QC Sonia Khillan 27SEP2023
Name Title Signature/Date

NF-light® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.517
C	1.59
D	5.17
E	15.1
F	50.4
G	149
H	431

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	7.66
C	24.2
D	78.8
E	237
F	880
G	2507
H	6673