

Product Number:	104080
Lot Number:	503881
Expiration:	9-May-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	312903	312903A
Detector Reagent	103512	315203	315203A
SBG Reagent	103514	315204	315204A
Sample Diluent	103516	315107	315107A
Calibrators	103517	313004	313004A
Control 0	104079	313104	313104A
Control 1	104078	313104	313104A
Control 2	104077	313104	313104A
RGP Reagent	103159	N/A <sup>1</sup>	N/A <sup>1</sup>

<sup>1</sup> 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	8.45	6.76 - 10.1
Control 1	313104	149	106 - 159
Control 2	313104	415	313 - 470
Panel 1	310007	7.57	4.60 - 9.38
Panel 2	310007	4.28	2.44 - 6.04
Panel 3	306904	98.0	74.0 - 111
Panel 4	306904	526	359 - 597

**GFAP®**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 0	313104	75.1	49.3 - 86.8
Control 1	313104	432	333 - 540
Control 2	313104	1924	1671 - 2506
Panel 1	310007	68.9	45.6 - 86.0
Panel 2	310007	63.0	40.4 - 73.7
Panel 3	306904	1080	795 - 1440
Panel 4	306904	7098	6199 - 9299

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

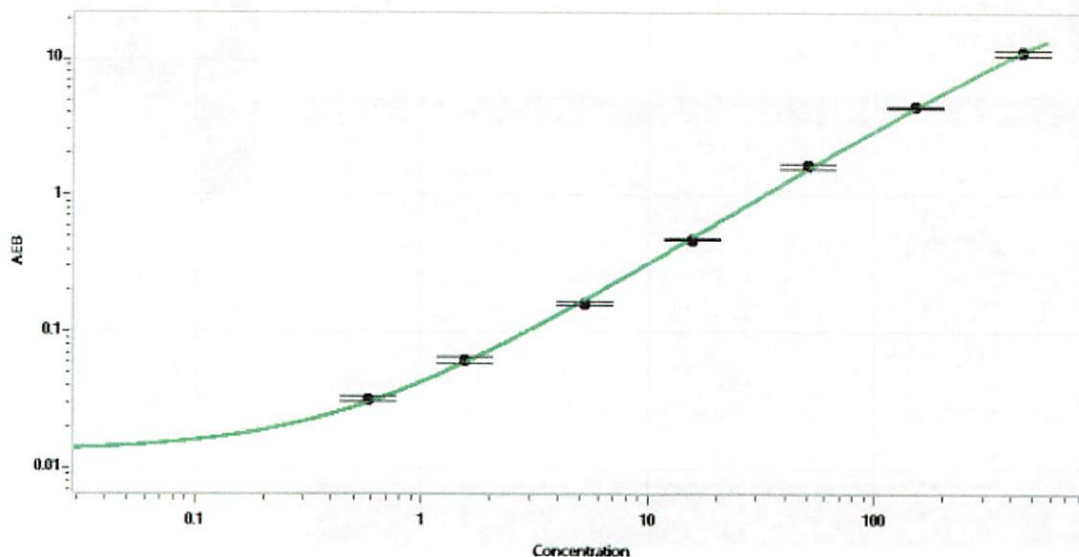
Review/Approval

Sonia Khillan  
Name

QC Manager  
Title

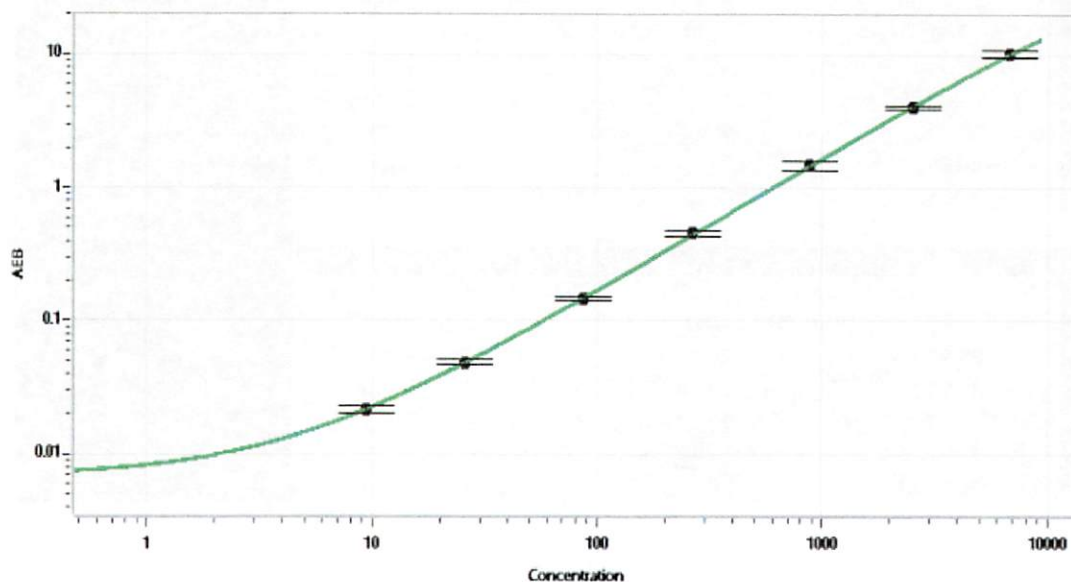
Sonia Khillan 13 JUN 2022  
Signature/Date

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