

Product Number:	103520
Lot Number:	503999
Expiration:	2-Sep-2024
Platform(s):	HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	103510	328413
Detector Reagent	103512	328510
SBG Reagent	103514	330402
Sample Diluent	103516	328512
Calibrators	103517	321906
Control 1	103518	321907
Control 2	103519	321907
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	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 1	321907	6.11	4.87 - 7.30
Control 2	321907	536	438 - 657
Panel 1	310007	6.27	4.40 - 11.0
Panel 2	323304	5.53	4.46 - 8.31
Panel 3	306904	76.9	74.0 - 111
Panel 4	306904	399	359 - 597

GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	321907	108	61.0 - 117
Control 2	321907	9158	6107 - 10721
Panel 1	310007	74.2	60.3 - 103
Panel 2	323304	60.6	34.3 - 82.1
Panel 3	306904	1204	795 - 1440
Panel 4	306904	8100	6199 - 9299

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

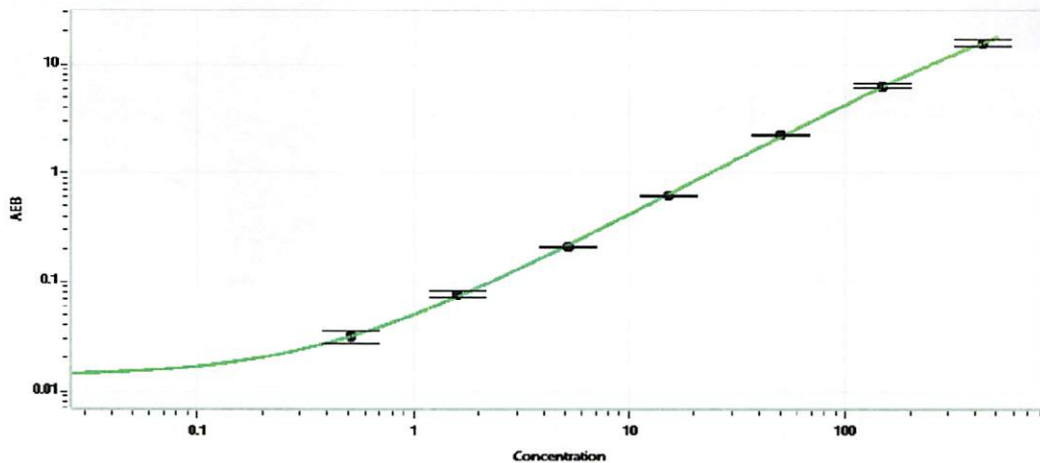
Review/Approval

Sonia Khillan
Name

QC
Title

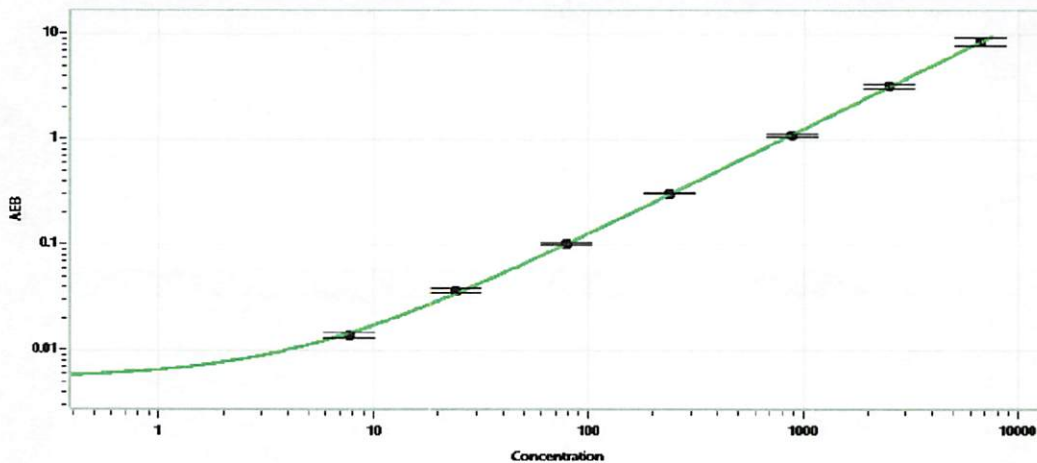
Sonia Rella 05 DEC 2023
Signature/Date

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	0.00
B	7.66
C	24.2
D	78.8
E	237
F	880
G	2507
H	6673