

Product Number:	103520
Lot Number:	503858
Expiration:	21-Jun-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	103510	317211	317211A
Detector Reagent	103512	321313	N/A
SBG Reagent	103514	320905	N/A
Sample Diluent	103516	317307	317307A
Calibrators	103517	320202	N/A
Control 1	103518	320203	N/A
Control 2	103519	320203	N/A
SBG Diluent (SR-X)†	103158	N/A ¹	N/A ¹
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 1	320203	7.22	5.56 - 8.34
Control 2	320203	539	439 - 692
Panel 1	310007	7.78	4.60 - 9.38
Panel 2	323304	6.68	4.60 - 7.45
Panel 3	306904	78.1	74.0 - 111
Panel 4	306904	405	359 - 597

GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	320203	104	82.4 - 137
Control 2	320203	11143	8812 - 13219
Panel 1	310007	79.1	45.6 - 86.0
Panel 2	323304	63.6	31.8 - 77.7
Panel 3	306904	1109	795 - 1440
Panel 4	306904	7883	6199 - 9299

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

Review/Approval

Sonia Khillan

Name

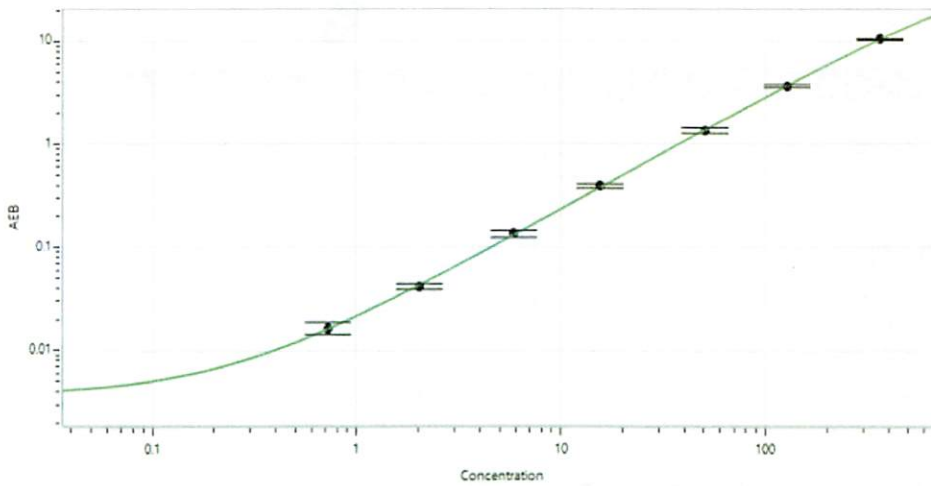
QC

Title

Sonia Rella 07DEC2023

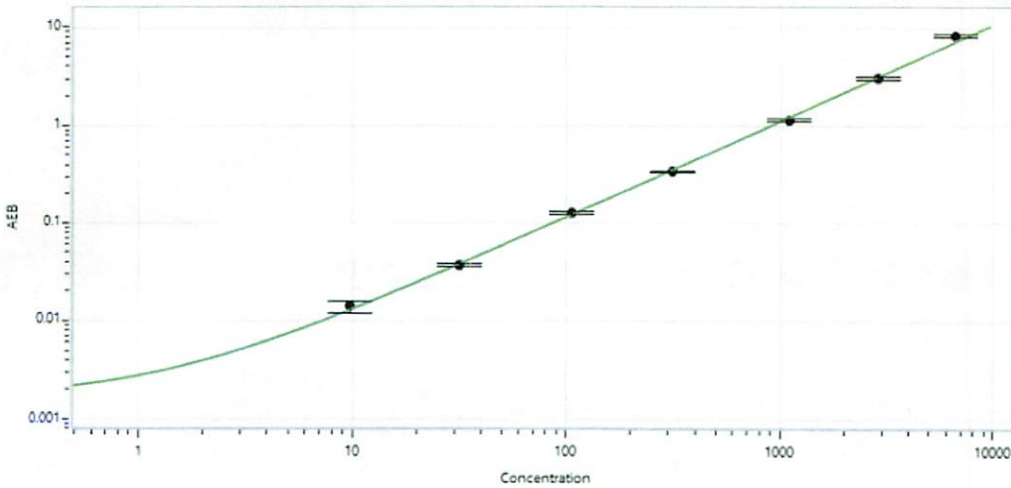
Signature/Date

NF-light® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.677
C	1.84
D	5.70
E	15.6
F	52.2
G	139
H	394

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	10.0
C	30.5
D	108
E	316
F	1115
G	3008
H	7350