

Product Number:	104080
Lot Number:	503797
Expiration:	11-Nov-2023
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103510	231507
Detector Reagent	103512	302703
SBG Reagent	103514	232007
Sample Diluent	103516	232008
Calibrators	103517	221303 / 224903
Control 0	104079	304110
Control 1	104078	304110
Control 2	104077	304110
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 0	304110	3.61	2.24 - 4.83
Control 1	304110	74.7	56.3 - 84.4
Control 2	304110	221	170 - 255
Panel 1	207003	13.3	10.7 - 16.1
Panel 2	207003	6.09	4.77 - 7.68
Panel 3	207003	134	101 - 151
Panel 4	207003	349	273 - 416

GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 0	304110	67.2	52.8 - 80.1
Control 1	304110	432	337 - 505
Control 2	304110	2080	1677 - 2515
Panel 1	207003	56.6	49.4 - 83.9
Panel 2	207003	65.2	52.9 - 104
Panel 3	207003	980	684 - 1104
Panel 4	207003	5657	4259 - 7215

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

Review/Approval

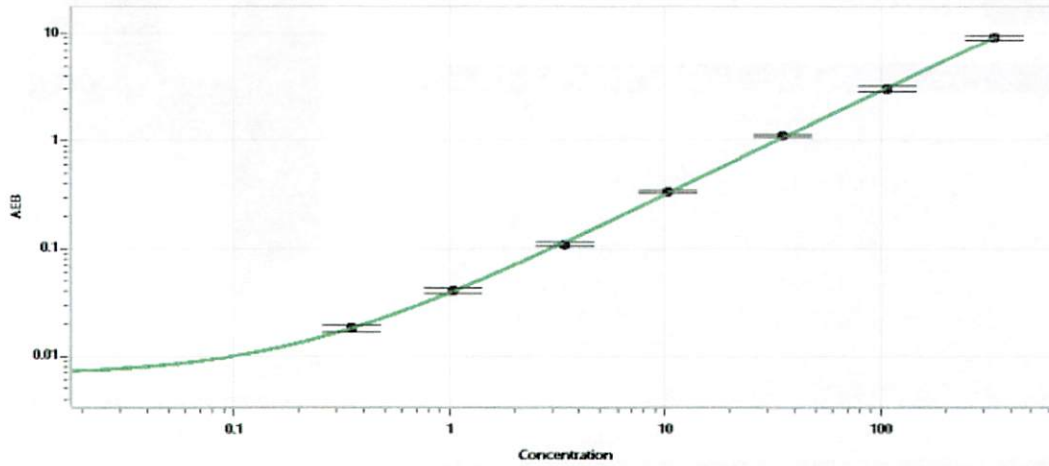
Sonia Khilkin
Name

QC

Title

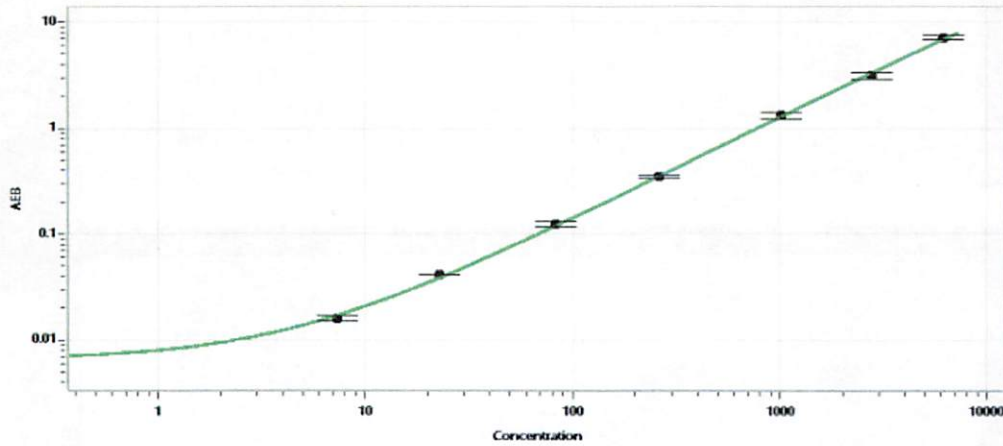
Sonia Khilkin 03MAR2023
Signature/Date

NF-light® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.352
C	1.04
D	3.42
E	10.3
F	35.3
G	107
H	335

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	7.36
C	22.9
D	82.5
E	261
F	1014
G	2766
H	6154