

Product Number:	103520
Lot Number:	503812
Expiration:	17-Feb-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	305205
Detector Reagent	103512	305315
SBG Reagent	103514	307304
Sample Diluent	103516	305308
Calibrators	103517	305807
Control 1	103518	305808
Control 2	103519	305808
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	305808	7.45	4.45 - 8.21
Control 2	305808	507	403 - 604
Panel 1	310007	8.60	4.21 - 9.64
Panel 2	310007	5.46	2.12 - 6.42
Panel 3	306904	93.9	73.9 - 110.8
Panel 4	306904	412	362 - 594

GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	305808	129	79.2 - 144
Control 2	305808	11758	9066 - 13598
Panel 1	310007	56.8	43.2 - 85.0
Panel 2	310007	50.9	41.1 - 69.8
Panel 3	306904	1320	742 - 1457
Panel 4	306904	7908	6356 - 9535

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

Review/Approval

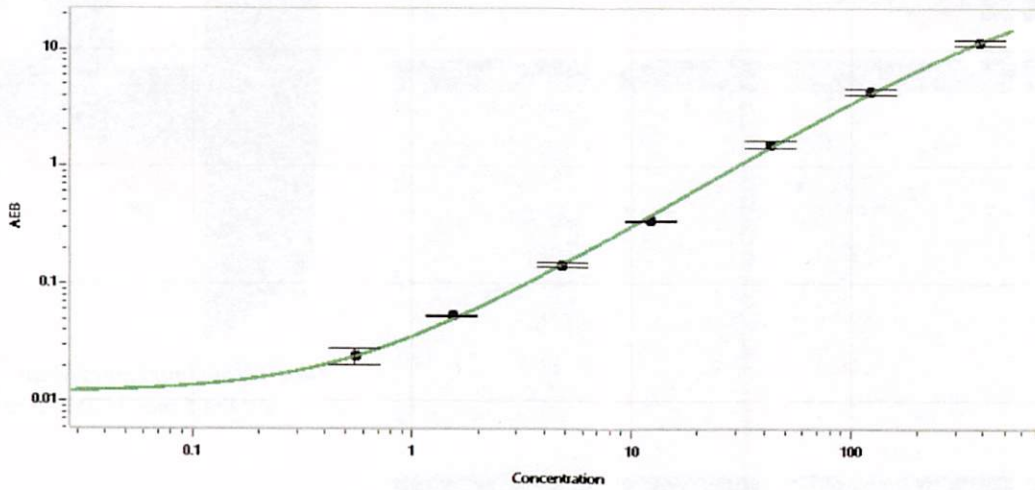
Sonia Khillan
Name

QC

Title

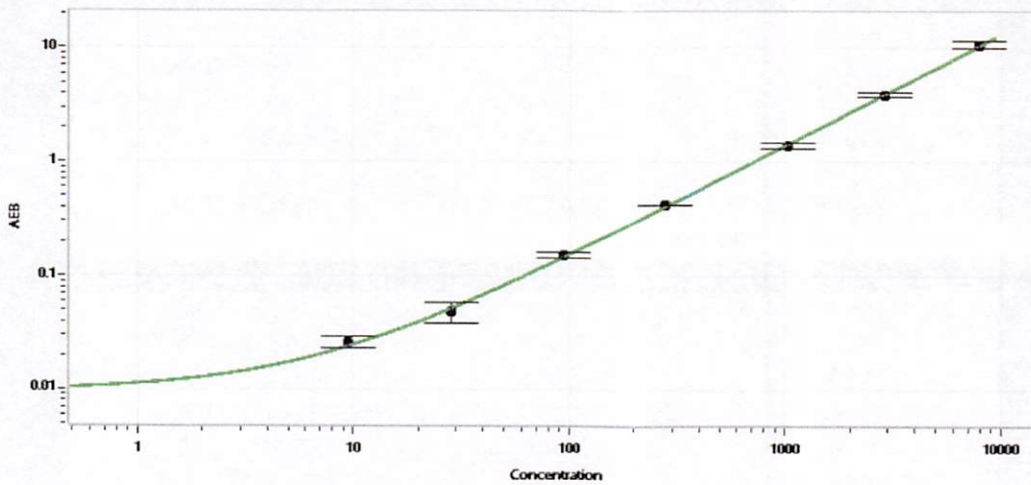
Sonia Khillan 27 APR 2023
Signature/Date

NF-light® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.552
C	1.53
D	4.81
E	12.2
F	42.7
G	122
H	387

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	9.54
C	28.5
D	93.9
E	279
F	1029
G	2898
H	7883