

Product Number:	103520
Lot Number:	503445
Expiration:	18-Feb-2023
Platform(s):	HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103510	204906
Detector Reagent	103512	204907
SBG Reagent	103514	210206
Sample Diluent	103516	205002
Calibrators	103517	221303/224903
Control 1	103518	221505
Control 2	103519	221505
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹



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

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	221505	4.51	3.07-5.50
Control 2	221505	418	336-505
Panel 1	207003	13.9	10.7-16.1
Panel 2	207003	6.69	4.77-7.68
Panel 3	207003	125	101-151
Panel 4	207003	328	273-416

GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	221505	96.4	74.7-112
Control 2	221505	10677	8118-12177
Panel 1	207003	62.4	49.4-83.9
Panel 2	207003	75.2	52.9-104
Panel 3	207003	920	684-1104
Panel 4	207003	5808	4259-7215

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

Review/Approval

Sonia Khillan

Name

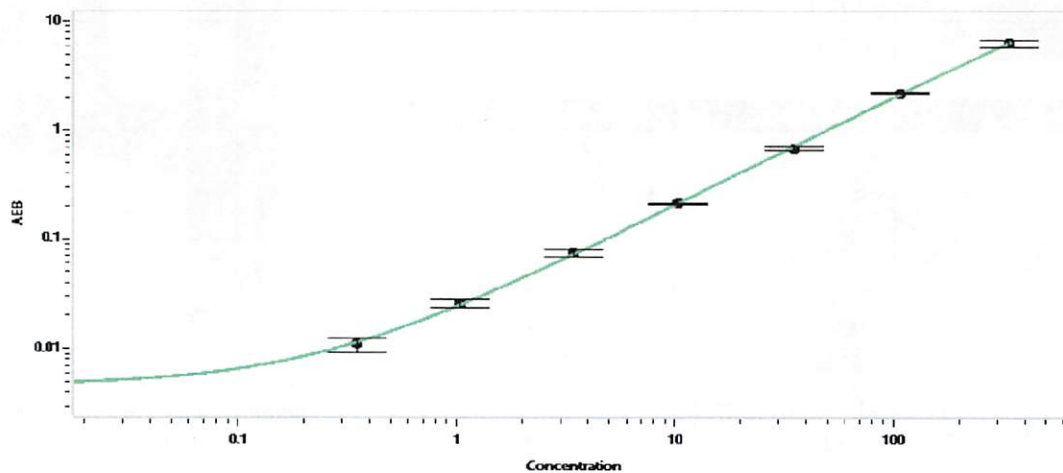
Ac Manager

Title

Sonia Khillan / 20SEP2022

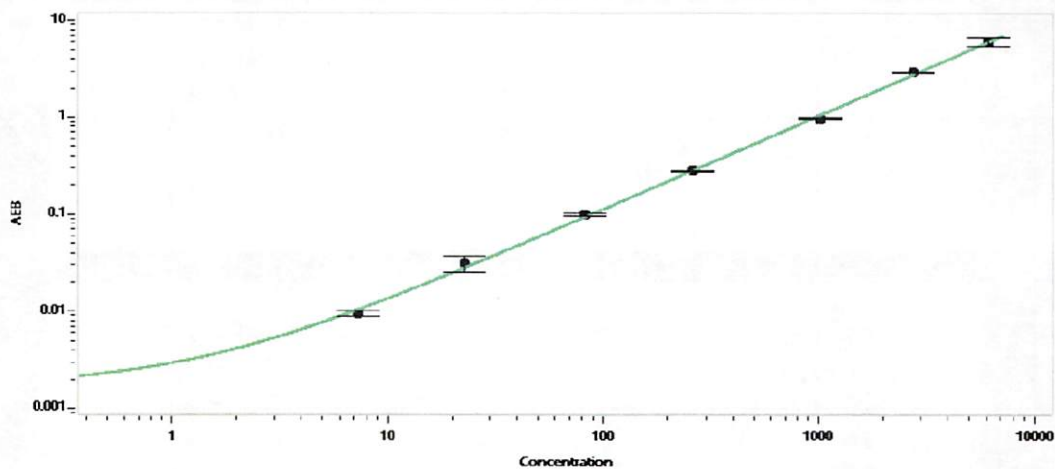
Signature/Date

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