

Product Number:	103520
Lot Number:	503514
Expiration:	13-Jul-2023
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	219407
Detector Reagent	103512	219904
SBG Reagent	103514	226501
Sample Diluent	103516	219906
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

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	3.60	2.87-5.47
Control 2	221505	397	335-502
Panel 1	207003	13.8	10.7-16.1
Panel 2	207003	6.42	4.77-7.68
Panel 3	207003	131	101-151
Panel 4	207003	386	273-416

GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	221505	94.3	73.6-111
Control 2	221505	9535	8061-12092
Panel 1	207003	57.4	49.4-83.9
Panel 2	207003	70.1	52.9-104
Panel 3	207003	945	684-1104
Panel 4	207003	5933	4259-7215

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

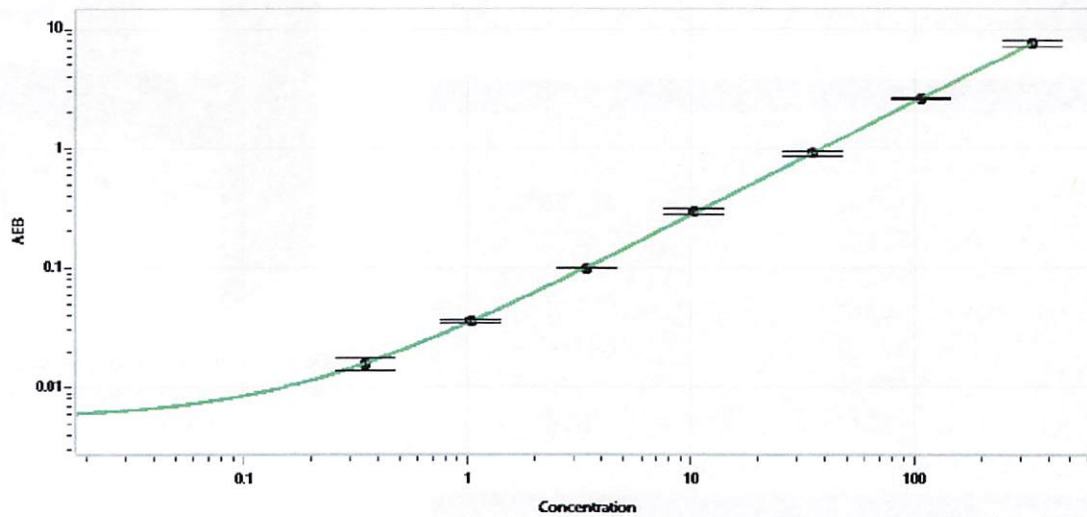
Review/Approval

Sonia Khillan
Name

QC Manager
Title

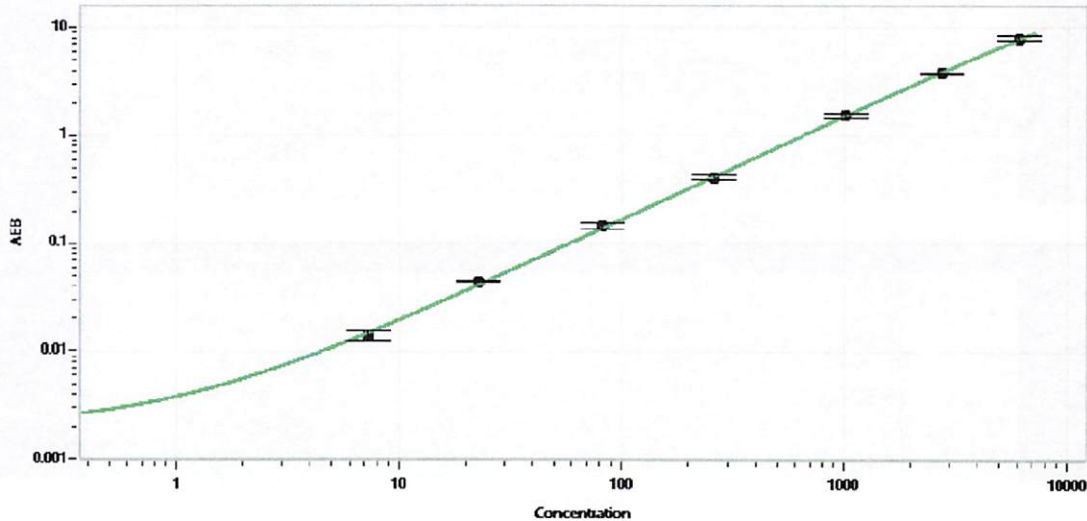
Sonia Khillan / 18 Oct 2022
Signature/Date

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