

Product Number:	103520
Lot Number:	503754
Expiration:	14-Nov-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	233904
Detector Reagent	103512	234006
SBG Reagent	103514	234007
Sample Diluent	103516	234010
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	3.92	2.87 - 5.47
Control 2	221505	382	335 - 502
Panel 1	207003	15.6	10.7 - 16.1
Panel 2	207003	7.19	4.77 - 7.68
Panel 3	207003	138	101 - 151
Panel 4	207003	408	273 - 416

GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	221505	95.5	73.6 - 111
Control 2	221505	9952	8061 - 12092
Panel 1	207003	70.9	49.4 - 83.9
Panel 2	207003	82.5	52.9 - 104
Panel 3	207003	992	684 - 1104
Panel 4	207003	6370	4259 - 7215

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

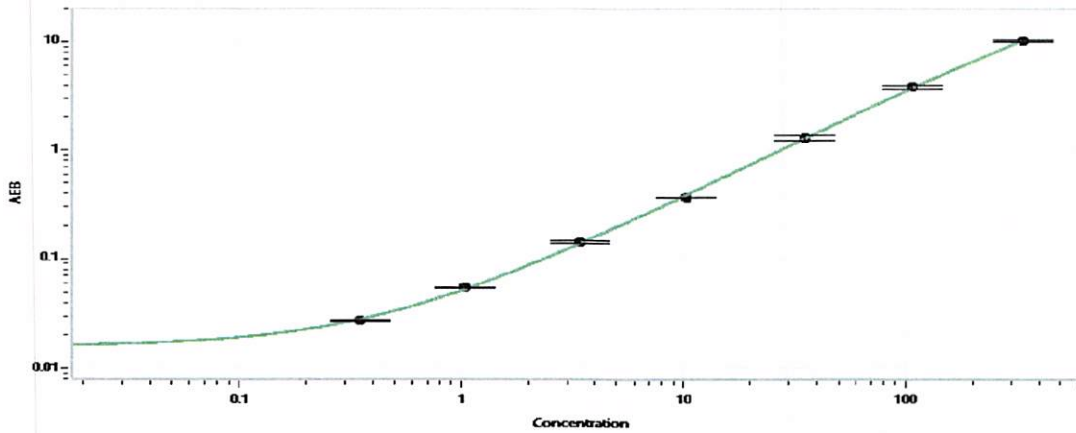
Review/Approval

Sonia Khillan
Name

QC
Title

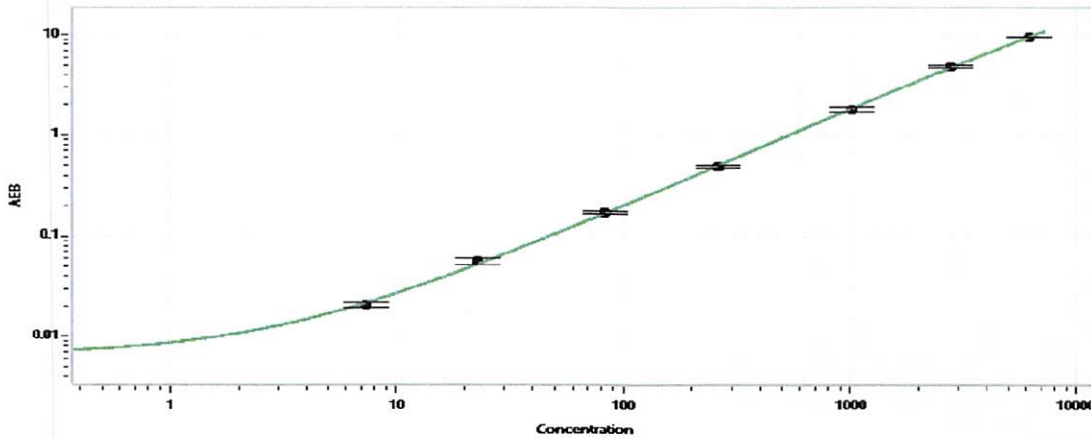
Sonia Khillan 16DEC2022
Signature/Date

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