

Product Number:	104080
Lot Number:	503518
Expiration:	20-Apr-2023
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103510	204906
Detector Reagent	103512	204907
SBG Reagent	103514	205001
Sample Diluent	103516	205002
Calibrators	103517	214507
Control 0	104079	214601
Control 1	104078	214601
Control 2	104077	214601
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	214601	8.00	5.20-9.09
Control 1	214601	78.5	61.0-91.5
Control 2	214601	231	182-272
Panel 1	207003	14.1	10.7-16.1
Panel 2	207003	6.34	4.77-7.68
Panel 3	207003	130	101-151
Panel 4	207003	342	273-416

GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 0	214601	67.4	50.3-79.7
Control 1	214601	395	305-458
Control 2	214601	1906	1580-2369
Panel 1	207003	70.2	49.4-83.9
Panel 2	207003	69.1	52.9-104
Panel 3	207003	1030	684-1104
Panel 4	207003	5942	4259-7215

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

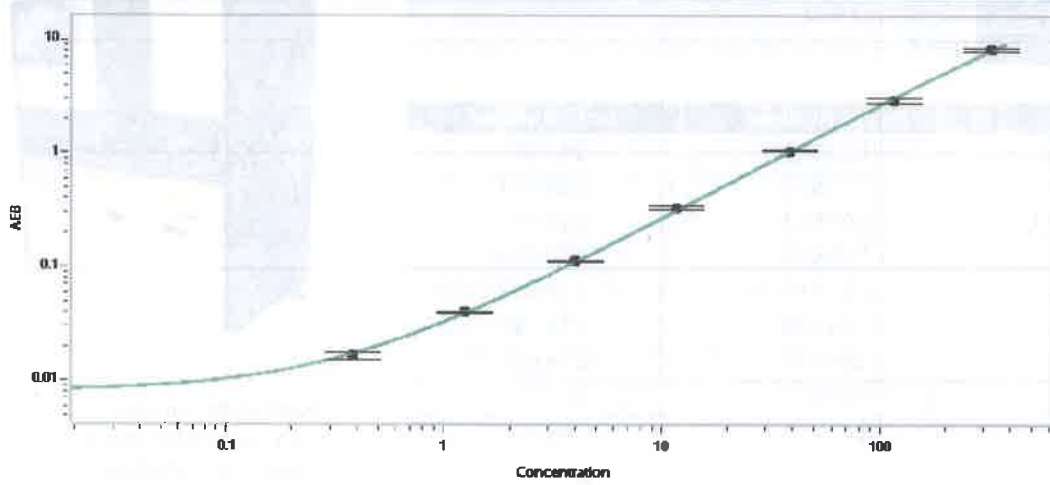
Review/Approval

Sonia Khillan
Name

QC Manager
Title

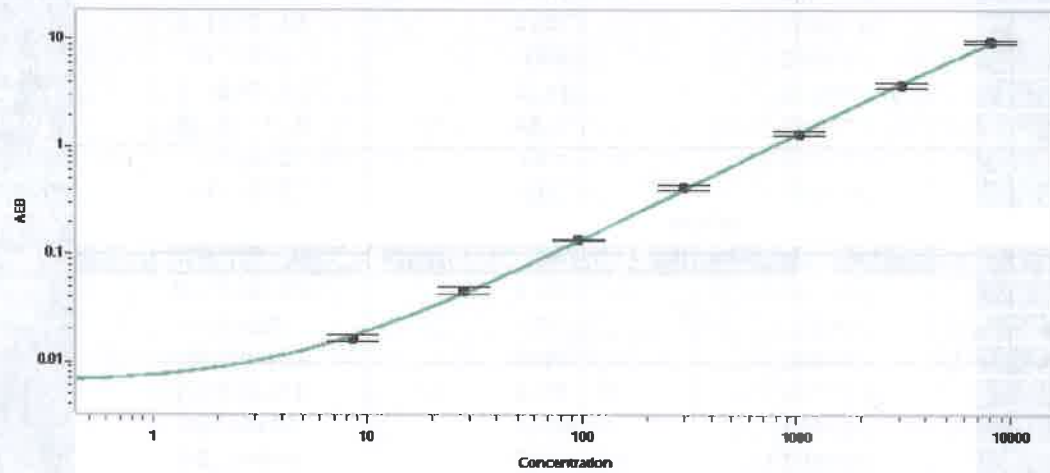
Sonia Khillan - 07JUL2022
Signature/Date

NF-light® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.383
C	1.24
D	4.02
E	11.7
F	38.5
G	116
H	323

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	8.60
C	28.3
D	95.6
E	295
F	1031
G	3072
H	8008