

Product Number:	103520
Lot Number:	502553
Expiration:	2-Mar-2021
Platform(s):	HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	103510	006203
Detector Reagent	103512	006515
SBG Reagent	103514	006201
Sample Diluent	103516	027411
Calibrators	103517	008701
Control 1	103518	008702
Control 2	103519	008702
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is 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).



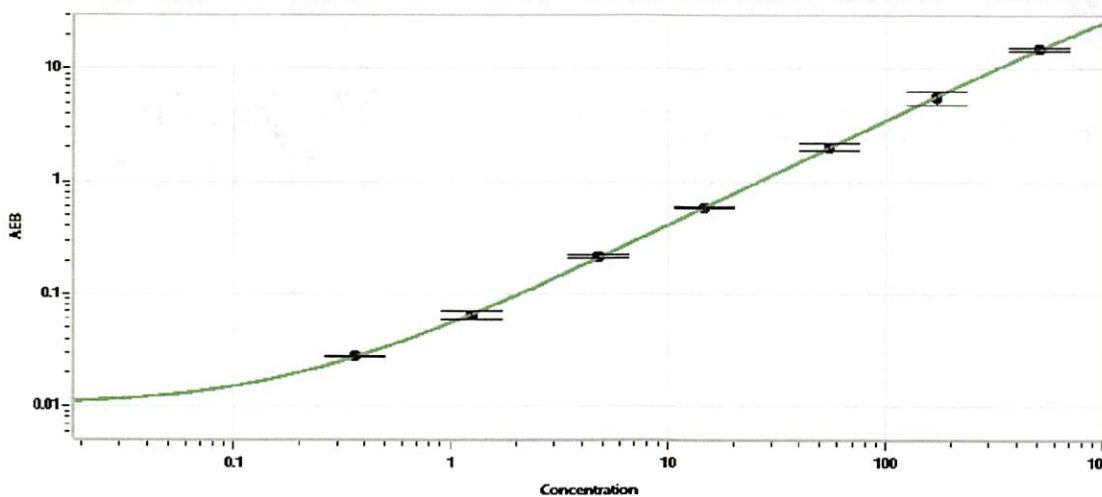
Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials			
NF-light®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	008702	4.49	3.23-7.95
Control 2	008702	581	495-875
Panel 1	929401	3.07	2.56-4.94
Panel 2	929402	6.12	5.09-11.1
Panel 3	929403	98.3	77.5-116
Panel 4	929404	473	375-562
GFAP®			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	008702	74.9	73.9-111
Control 2	008702	10674	9108-14850
Panel 1	929401	45.2	38.9-66.8
Panel 2	929402	44.0	39.3-74.2
Panel 3	929403	1304	1028-1542
Panel 4	929404	4599	3500-5314

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

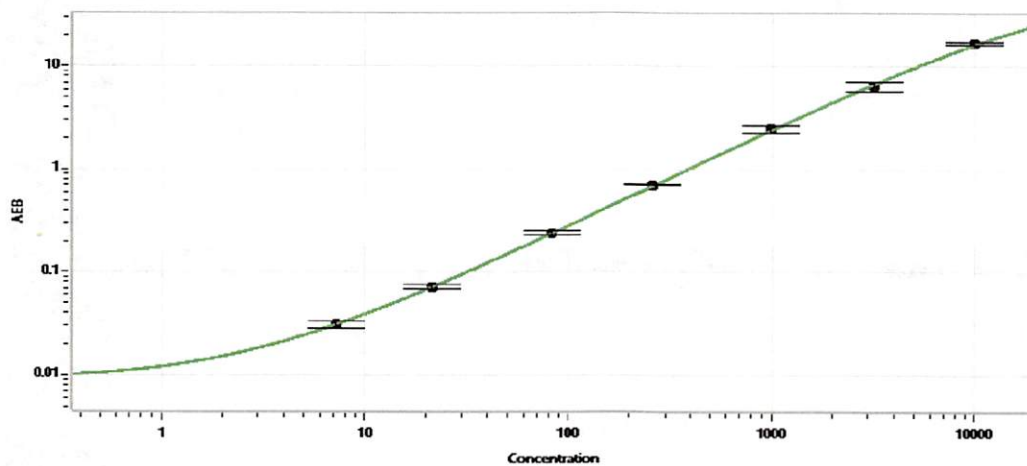
Review/Approval	Shivani Goel	QC Manager	Shivani Goel 20 Oct 2020
	Name	Title	Signature/Date

NF-light® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.364
C	1.25
D	4.75
E	14.5
F	54.6
G	171
H	509

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	7.17
C	21.6
D	83.7
E	260
F	990
G	3175
H	9950