

Product Number:	103520
Lot Number:	502816
Expiration:	13-Mar-2022
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	107201
Detector Reagent	103512	107712
SBG Reagent	103514	107301
Sample Diluent	103516	107302
Calibrators	103517	107402
Control 1	103518	107403
Control 2	103519	107403
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).

Release Materials

NF-light®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	107403	7.28	5.64-8.45
Control 2	107403	672	530-795
Panel 1	929401	3.45	2.63-4.73
Panel 2	929402	6.74	4.56-10.5
Panel 3	929403	81.9	74.3-112
Panel 4	929404	366	325-554

GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	107403	116	85.2-129
Control 2	107403	10625	8745-13118
Panel 1	929401	61.1	39.8-64.9
Panel 2	929402	61.4	34.1-71.7
Panel 3	929403	1173	1007-1511
Panel 4	929404	4067	3421-5189

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

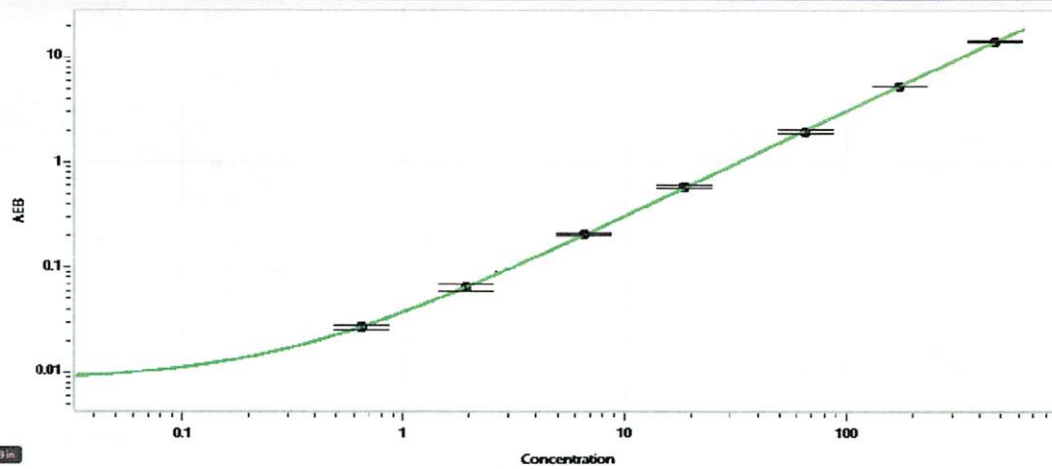
Review/Approval

Junpil Park
Name

QC Supervisor
Title

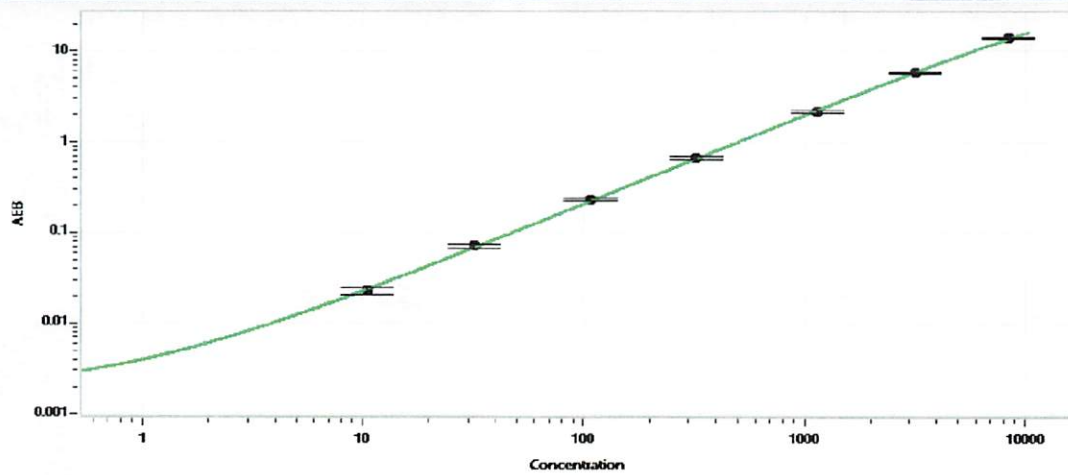

Signature/Date
17 May 2021

NF-light® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.646
C	1.93
D	6.54
E	18.7
F	65.1
G	175
H	473

GFAP® Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	10.5
C	31.9
D	108
E	322
F	1147
G	3154
H	8418