

Product Number:	103520
Lot Number:	503227
Expiration:	3-Sep-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	132404
Detector Reagent	103512	132405
SBG Reagent	103514	132507
Sample Diluent	103516	132407
Calibrators	103517	133406
Control 1	103518	133407
Control 2	103519	133407
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	133407	3.91	2.81 - 4.97
Control 2	133407	424	308 - 483
Panel 1	128902	9.94	8.62 - 14.0
Panel 2	128902	11.5	9.38 - 14.6
Panel 3	128902	115	99.4 - 149
Panel 4	128902	292	257 - 422

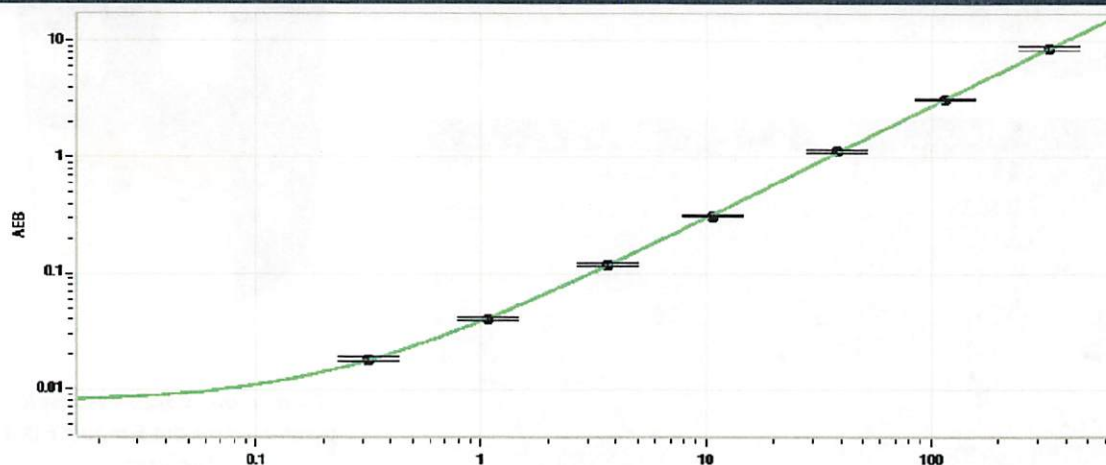
GFAP®

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	133407	108	68.6 - 129
Control 2	133407	11480	9072 - 13608
Panel 1	128902	78.9	46.7 - 95.3
Panel 2	128902	63.7	40.1 - 89.1
Panel 3	128902	5305	3360 - 5866
Panel 4	128902	21223	16761 - 25142

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

Review/Approval	Junpil Park	QC Supervisor	 14 Dec 2021
	Name	Title	Signature/Date

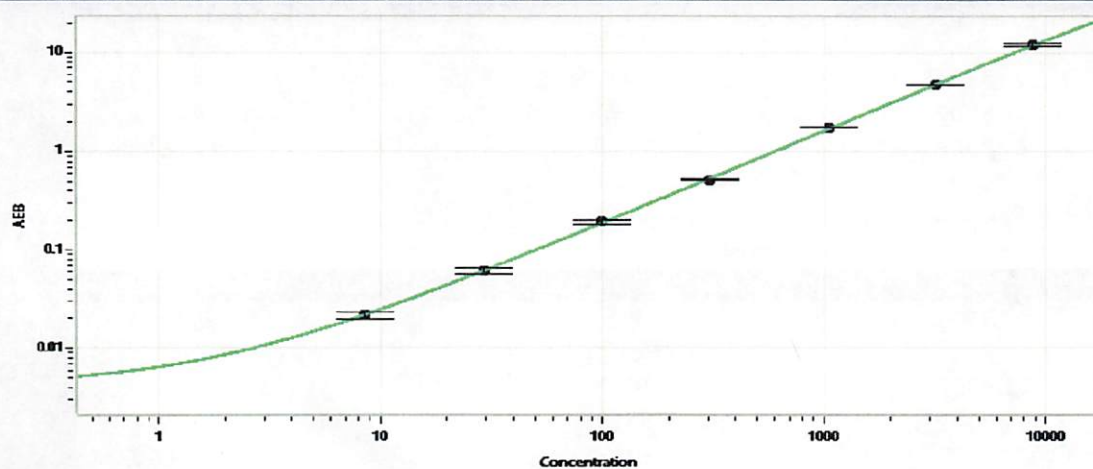
NF-light® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.318
C	1.07
D	3.65
E	10.7
F	38.3
G	116
H	336

GFAP® Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	8.47
C	29.2
D	99.9
E	307
F	1060
G	3170
H	8733