

Product Number:	103520
Lot Number:	504401
Expiration:	21-Feb-2026
Platform(s):	SR-X and HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103510	427403
Detector Reagent	103512	428113
SBG Reagent	103514	427704
Sample Diluent Reagent	103516	427601
Reference Calibrators	103517	428103
Control 1	103518	428104
Control 2	103519	428104
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

† Specific to use on the SR-X platform - refer to Neuro 2-Plex B SR-X Kit Instructions (KI-0062).

Release Materials*

NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	428104	5.78	3.94 - 6.63
Control 2	428104	497	369 - 553

GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	428104	86.6	65.5 - 98.2
Control 2	428104	8813	6892 - 10338

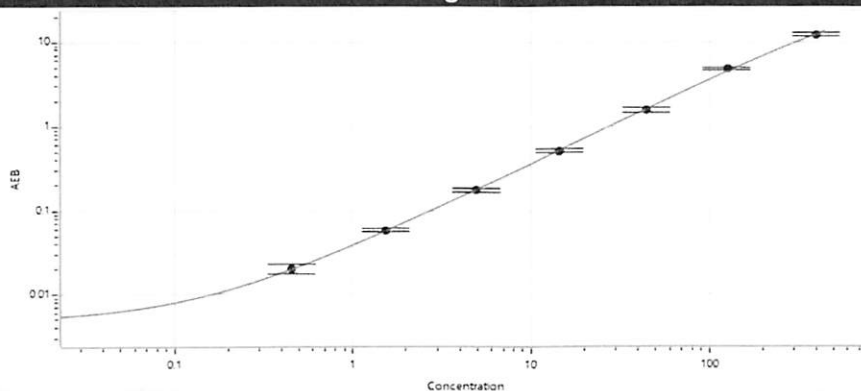
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Review/Approval

Dhvani Patel Doc Control I, QA DP 30 Oct 2024

Name Title Signature/Date

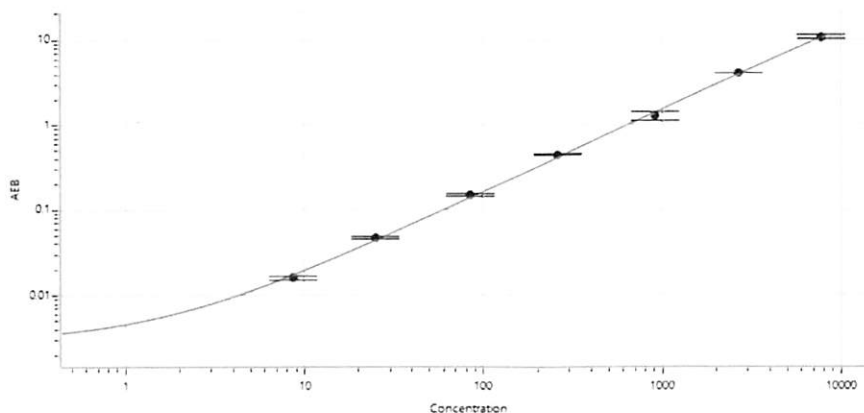
NF-Light® Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.457
C	1.54
D	4.93
E	14.4
F	44.4
G	127
H	401

GFAP® Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	8.61
C	25.0
D	84.9
E	261
F	910
G	2693
H	7777