

Product Number:	103520
Lot Number:	504653
Expiration:	13-Dec-2026
Platform(s):	SR-X and HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	103510	518402
Detector Reagent	103512	518403
SBG Reagent	103514	518306
Sample Diluent Reagent	103516	518203
Reference Calibrators	103517	518804
Control 1	103518	518805
Control 2	103519	518805
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).



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

Release Materials*			
NF-Light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	518805	5.68	4.25 - 6.85
Control 2	518805	496	371 - 557

GFAP®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	518805	84.0	62.5 - 93.8
Control 2	518805	8946	6784 - 10177

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

Review/Approval

Shwami Patel

Name

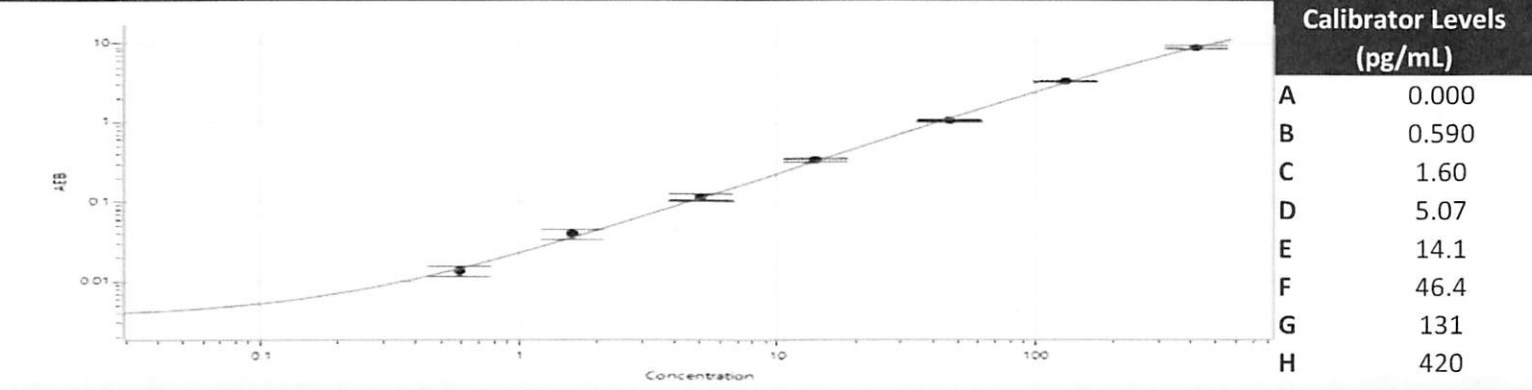
DOC Control/II, QA

Title

SPR / 08 AUG 2025

Signature/Date

NF-Light® Calibrator Curve



GFAP® Calibrator Curve

