

Product Number:	104670
Lot Number:	504629
Expiration:	23-Nov-2026
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	104652	515303	N/A
Detector Reagent	104654	515401	N/A
SBG Reagent	104656	515402	N/A
Plasma Sample Diluent	104722	514707	N/A
CSF Sample Diluent	104723	514314	N/A
Calibrator Diluent	104657	514312	N/A
Control Diluent	104721	514313	N/A
Calibrator Concentrate	104666	317341	N/A
Control 1 Concentrate	104668	317342	N/A
Control 2 Concentrate	104669	317343	317343A
RGP 2.0 Reagent	104321	N/A ¹	N/A ¹
Activation Buffer Reagent	104324	N/A ¹	N/A ¹



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

¹ Reagents are not Kit Lot Specific

Review/Approval

D. Patel
Name

Sr. Doc. Control - QA
Title

22 Jul 2025
Signature/Date

Release Materials*

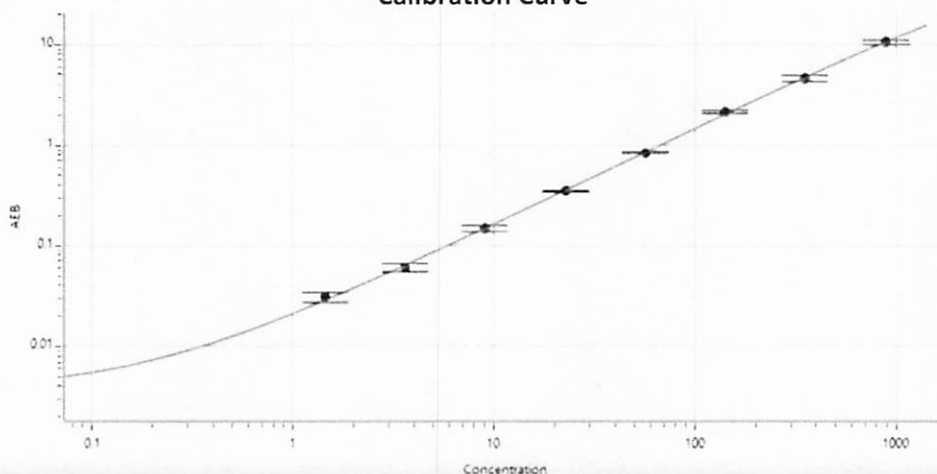
GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	317342	111	86.8 - 161
Control 2	317343A	1375	1049 - 1948

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

Calibrator Concentrate Reconstitution Volume (uL)	1144
Control 1 Concentrate Reconstitution Volume (uL)	1144
Control 2 Concentrate Reconstitution Volume (uL)	1144

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.45
C	3.62
D	9.03
E	22.6
F	56.6
G	142
H	354
I	885

Release Materials*

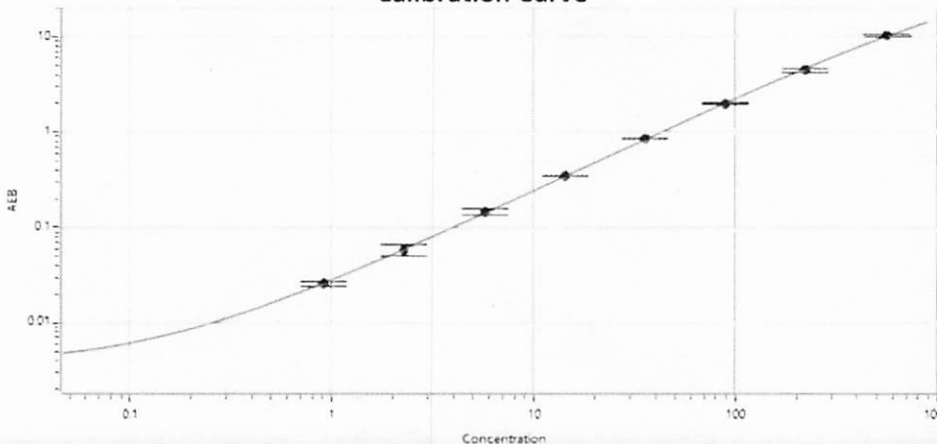
NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	317342	85.5	60.6 - 112
Control 2	317343A	1079	768 - 1427

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

Calibrator Concentrate Reconstitution Volume (uL)	1144
Control 1 Concentrate Reconstitution Volume (uL)	1144
Control 2 Concentrate Reconstitution Volume (uL)	1144

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.910
C	2.27
D	5.69
E	14.2
F	35.6
G	88.8
H	222
I	556