

Product Number:	104670
Lot Number:	504383
Expiration:	28-Feb-2026
Platform(s):	HD-X



Component	Part Number	Lot Number
Bead Reagent	104652	424805
Detector Reagent	104654	424901
SBG Reagent	104656	425603
Plasma Sample Diluent	104722	424310
CSF Sample Diluent	104723	424213
Calibrator Diluent	104657	424717
Control Diluent	104721	424718
Calibrator Concentrate	104666	402948
Control 1 Concentrate	104668	402949
Control 2 Concentrate	104669	402950
RGP 2.0 Reagent	104321	N/A ⁺
Activation Buffer Reagent	104324	N/A ⁺

⁺ Reagents are not Kit Lot Specific

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

Review/Approval

Dhruvi Patel

DOC Control I, QA

DP 08/11/2024

Name

Title

Signature/Date

Release Materials*

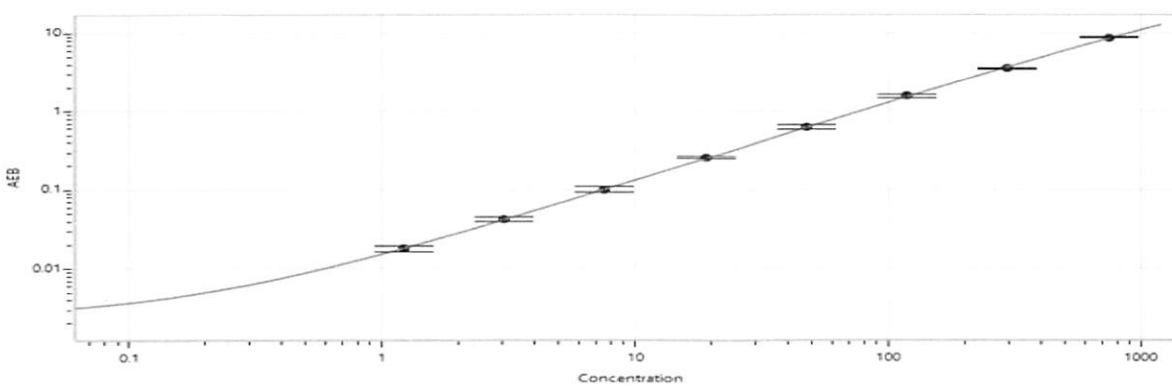
GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	402949	98.3	73.5 - 137
Control 2	402950	1242	855 - 1587

* 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	1.22
C	3.04
D	7.58
E	19.1
F	47.5
G	118
H	297
I	742

Release Materials*

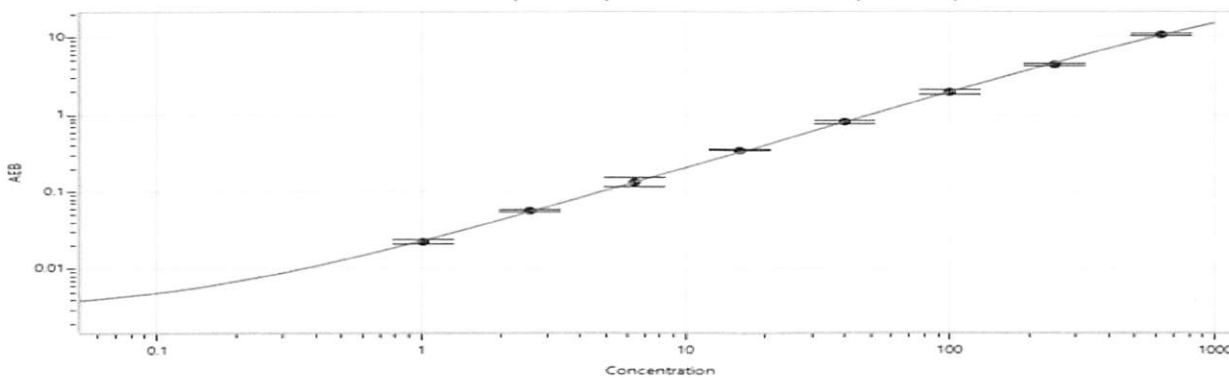
NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	402949	90.9	67.8 - 126
Control 2	402950	1179	798 - 1481

* 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	1.02
C	2.56
D	6.39
E	16.0
F	39.9
G	99.9
H	249
I	624