

<b>Product Number:</b>	104670
<b>Lot Number:</b>	504330
<b>Expiration:</b>	26-Jan-2026
<b>Platform(s):</b>	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	104652	421403
Detector Reagent	104654	421501
SBG Reagent	104656	421502
Plasma Sample Diluent	104722	421216
CSF Sample Diluent	104723	421117
Calibrator Diluent	104657	420808
Control Diluent	104721	420809
Calibrator Concentrate	104666	402913
Control 1 Concentrate	104668	402914
Control 2 Concentrate	104669	402915
RGP 2.0 Reagent	104321	N/A <sup>†</sup>
Activation Buffer Reagent	104324	N/A <sup>†</sup>

<sup>†</sup> Reagents are not Kit Lot Specific

Review/Approval

Dipesh Patel Doc Control - IT QA  
Name Title

[Signature] / 25-SEP-2024  
Signature/Date

**Release Materials\***

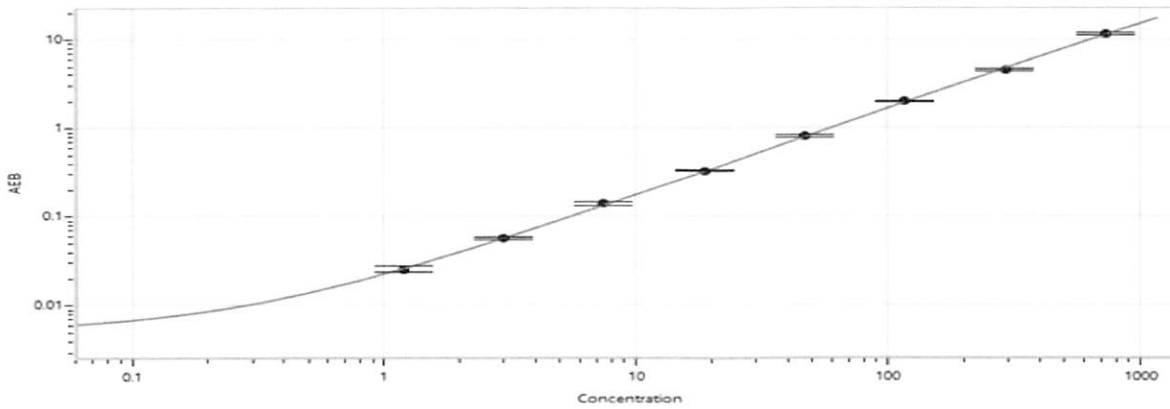
**GFAP®**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
<b>Control 1</b>	402914	101	70.7 - 131
<b>Control 2</b>	402915	1462	879 - 1633

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

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	1144
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	1144
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	1144

**Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	1.20
C	2.99
D	7.46
E	18.8
F	46.7
G	116
H	292
I	729

**Release Materials\***

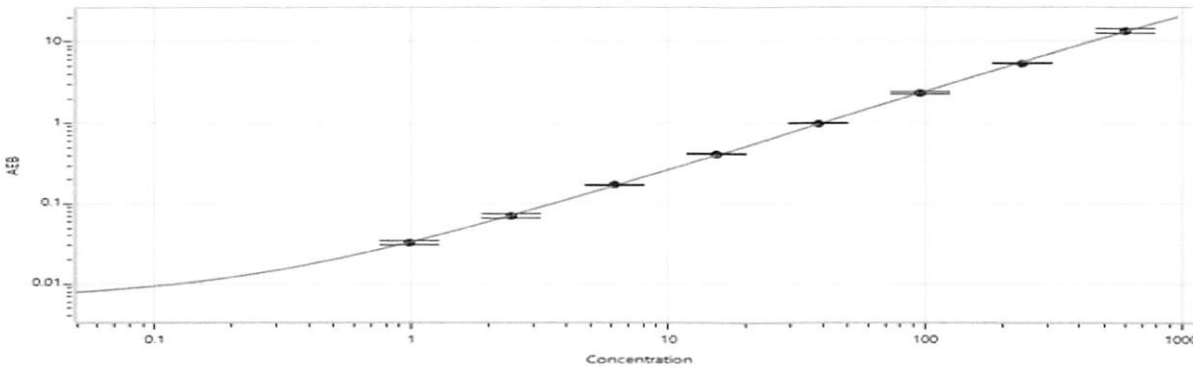
**NF-Light®**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
<b>Control 1</b>	402914	94.4	62.9 - 117
<b>Control 2</b>	402915	1142	804 - 1492

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

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	1144
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	1144
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	1144

**Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.983
C	2.46
D	6.14
E	15.4
F	38.3
G	96.0
H	239
I	600