

<b>Product Number:</b>	104670
<b>Lot Number:</b>	504092
<b>Expiration:</b>	7-Jan-2025
<b>Platform(s):</b>	HD-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	104652	334517	N/A
Detector Reagent	104654	334709	N/A
SBG Reagent	104656	334708	N/A
Plasma Sample Diluent	104722	329813	N/A
CSF Sample Diluent	104723	329626	N/A
Calibrator Diluent	104657	318815	318815A
Control Diluent	104721	318816	318816A
Calibrator Concentrate	104666	402913	N/A
Control 1 Concentrate	104668	402914	N/A
Control 2 Concentrate	104669	402915	N/A
RGP 2.0 Reagent	104321	N/A <sup>+</sup>	N/A <sup>+</sup>
Activation Buffer Reagent	104324	N/A <sup>+</sup>	N/A <sup>+</sup>

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

Review/Approval

Cody Carlson

QC Manager

Name

Title

*Cody Carlson* 02 APR 2024  
Signature/Date

**Release Materials\***

**GFAP®**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	402914	83.7	56.7 - 111
Control 2	402915	1208	872 - 1620
Panel 1	327811	63.8	48.3 - 89.7
Panel 2	327811	1624	1138 - 2113

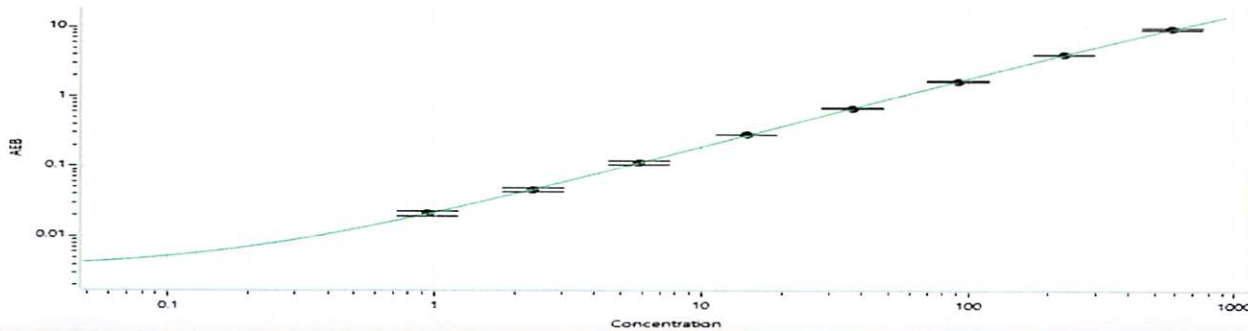
\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

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

**Calibrator Levels (pg/mL)**

A	0
B	0.944
C	2.36
D	5.88
E	14.8
F	36.8
G	91.7
H	230
I	575

**Calibration Curve**



**Release Materials\***

**NF-Light®**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	402914	78.6	48.2 - 106
Control 2	402915	1036	696 - 1292
Panel 1	327811	16.9	11.9 - 22.0
Panel 2	327811	464	310 - 575

\* 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

**Calibrator Levels (pg/mL)**

A	0
B	0.839
C	2.10
D	5.24
E	13.1
F	32.7
G	81.9
H	204
I	512

**Calibration Curve**

