

<b>Product Number:</b>	104822
<b>Lot Number:</b>	504356
<b>Expiration:</b>	12-Feb-2026
<b>Platform(s):</b>	HD-X



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

Component	Part Number	Lot Number	Control Number
Bead Reagent	104800	422902	N/A
Detector Reagent	104802	423302	N/A
SBG Reagent	104804	423303	N/A
Calibrator Diluent	104808	422622	N/A
Control Diluent	104809	422623	N/A
Plasma Sample Diluent	104805	422518	N/A
CSF Sample Diluent	104806	422521	N/A
Calibrator Concentrate	104818	317324	317324A
Control 1 Concentrate	104820	317325	317325A
Control 2 Concentrate	104821	317326	317326A
RGP 2.0 Reagent	104321	N/A <sup>1</sup>	N/A <sup>1</sup>
Activation Buffer Reagent	104324	N/A <sup>1</sup>	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific

Review/Approval

TAM LEE DOC CONTROL II, QA Man Zee 15 NOV 2024  
 Name Title Signature/Date

**Release Materials\***

**NF-Light®**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	317325A	19.4	14.9 - 27.9
Control 2	317326A	904	620 - 1151

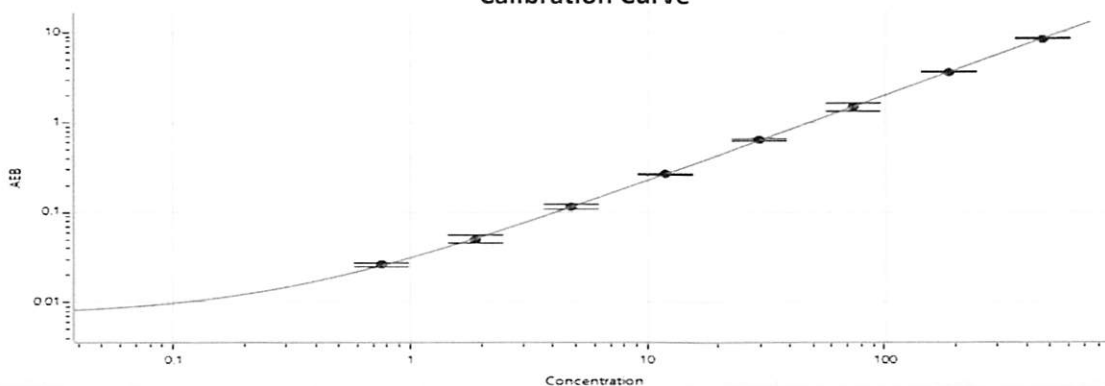
\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	1250
Control 1 Concentrate Reconstitution Volume (uL)	1250
Control 2 Concentrate Reconstitution Volume (uL)	1250

**Calibrator Levels (pg/mL)**

A	0.000
B	0.760
C	1.90
D	4.75
E	11.9
F	29.7
G	74.3
H	186
I	464

**Calibration Curve**



**Release Materials\***

**GFAP®**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	317325A	32.2	23.0 - 42.6
Control 2	317326A	1380	928 - 1723

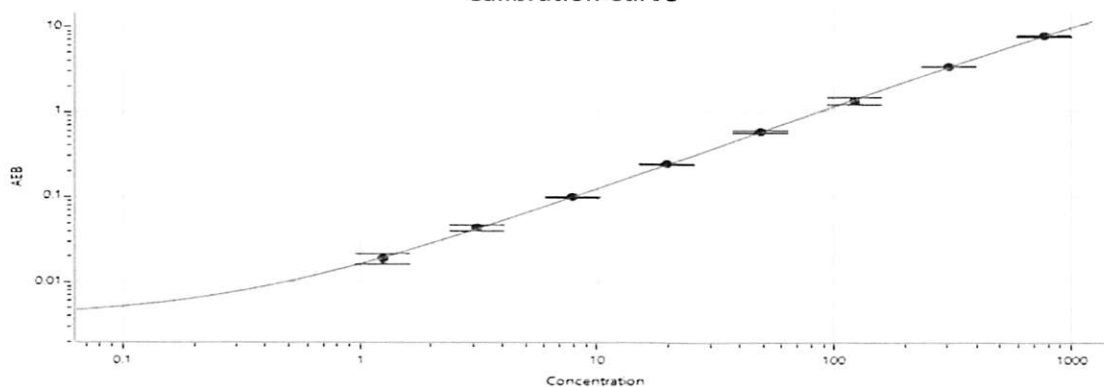
\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	1250
Control 1 Concentrate Reconstitution Volume (uL)	1250
Control 2 Concentrate Reconstitution Volume (uL)	1250

**Calibrator Levels (pg/mL)**

A	0.000
B	1.25
C	3.12
D	7.81
E	19.5
F	48.8
G	122
H	305
I	763

**Calibration Curve**



**Release Materials\***

**UCLH-1**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
<b>Control 1</b>	317325A	175	124 - 230
<b>Control 2</b>	317326A	6966	4737 - 8798

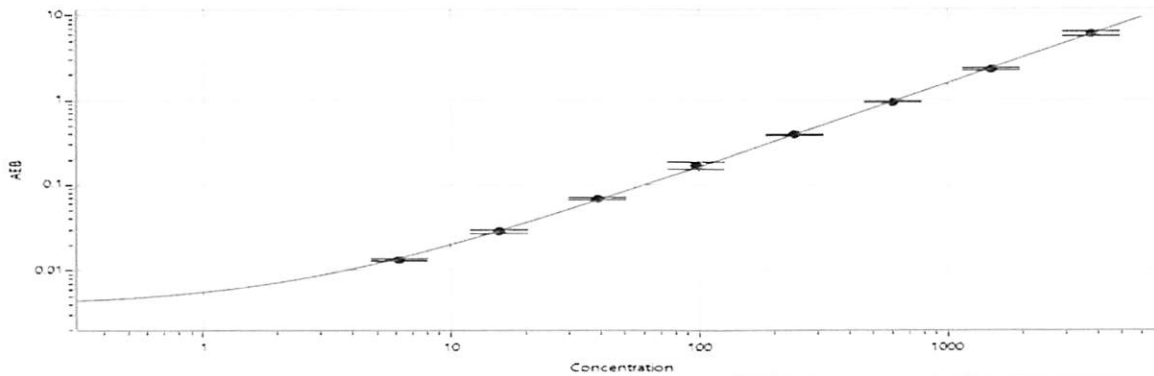
\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	1250
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	1250
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	1250

**Calibrator Levels (pg/mL)**

A	0.000
B	6.17
C	15.5
D	38.6
E	96.8
F	241
G	603
H	1507
I	3767

**Calibration Curve**



**Release Materials\***

**BD-Tau**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
<b>Control 1</b>	317325A	9.07	7.41 - 13.8
<b>Control 2</b>	317326A	434	284 - 528

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

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	1250
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	1250
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	1250

**Calibrator Levels (pg/mL)**

A	0.000
B	0.333
C	0.832
D	2.08
E	5.20
F	13.0
G	32.5
H	81.3
I	203

**Calibration Curve**

