

Product Number:	104822
Lot Number:	504408
Expiration:	24-Mar-2026
Platform(s):	HD-X



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

Component	Part Number	Lot Number	Control Number
Bead Reagent	104800	428406	N/A
Detector Reagent	104802	502901	502901A
SBG Reagent	104804	429006	N/A
Calibrator Diluent	104808	426816	N/A
Control Diluent	104809	426817	N/A
Plasma Sample Diluent	104805	428408	N/A
CSF Sample Diluent	104806	428509	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 ¹	N/A ¹
Activation Buffer Reagent	104324	N/A ¹	N/A ¹

¹ Reagents are not Kit Lot Specific

Review/Approval

Dhruvami Patel

Doc Control I, QA

DP 28 Feb 2025

Name

Title

Signature/Date

Release Materials*

NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	317325A	19.2	14.2 - 26.8
Control 2	317326A	948	585 - 1200

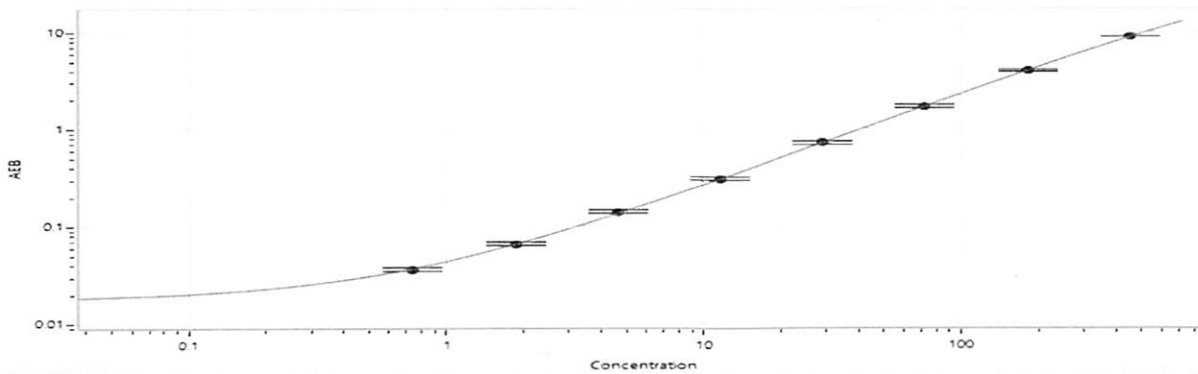
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

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	0.741
C	1.86
D	4.63
E	11.6
F	29.0
G	72.4
H	181
I	452

Calibration Curve



Release Materials*

GFAP®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	317325A	38.7	28.5 - 53.0
Control 2	317326A	1620	1105 - 2052

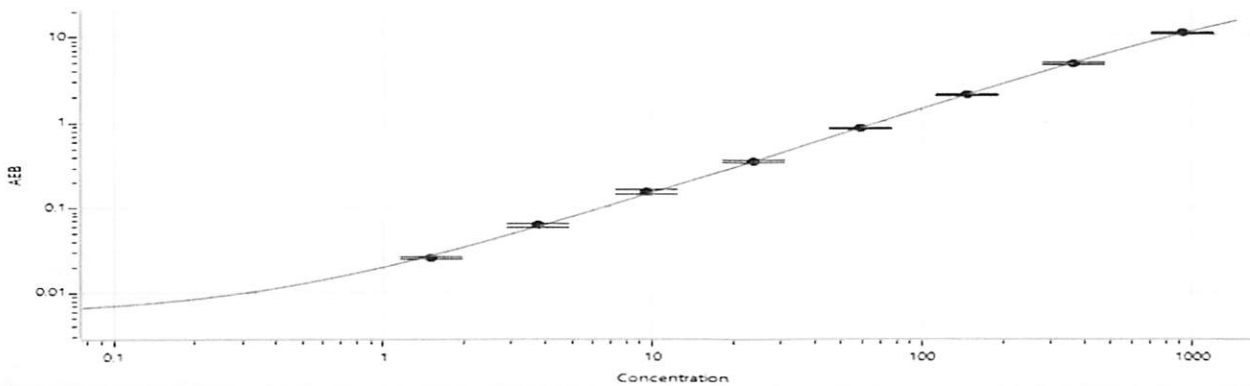
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

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	1.51
C	3.77
D	9.43
E	23.6
F	58.9
G	147
H	368
I	921

Calibration Curve



Release Materials*

UCLH-1

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	317325A	161	122 - 227
Control 2	317326A	6710	4991 - 9269

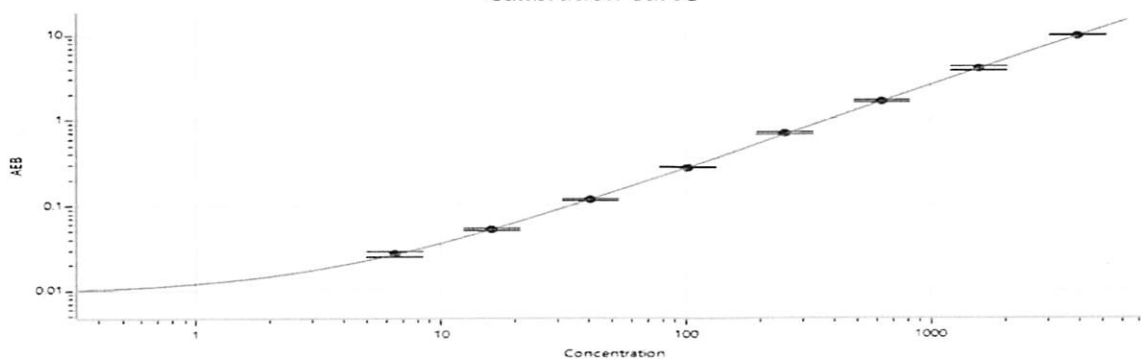
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

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	6.47
C	16.2
D	40.4
E	101
F	253
G	631
H	1578
I	3944

Calibration Curve



Release Materials*

BD-Tau

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	317325A	10.9	7.42 - 14.6
Control 2	317326A	438	301 - 559

* 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	0.331
C	0.827
D	2.07
E	5.17
F	13.0
G	32.3
H	80.8
I	202

Calibration Curve

