

<b>Product Number:</b>	104465
<b>Lot Number:</b>	504514
<b>Expiration:</b>	23-Mar-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	104448	426708	N/A
Detector Reagent	104450	426901	N/A
SBG Reagent	104452	503803	503803A
Calibrator Diluent Reagent	104453	426812	N/A
Plasma Sample Diluent Reagent	104647	426810	N/A
CSF Sample Diluent Reagent	104648	426809	N/A
Control Diluent Reagent	104649	426811	N/A
Calibrator Concentrate	104466	317328	317328A
Control 1 Concentrate	104468	317329	317329A
Control 2 Concentrate	104469	317330	317330A
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

*Dipesh Patel*

*Doc control TC*

*znt / 27 Feb 2025*

Name

Title

Signature/Date

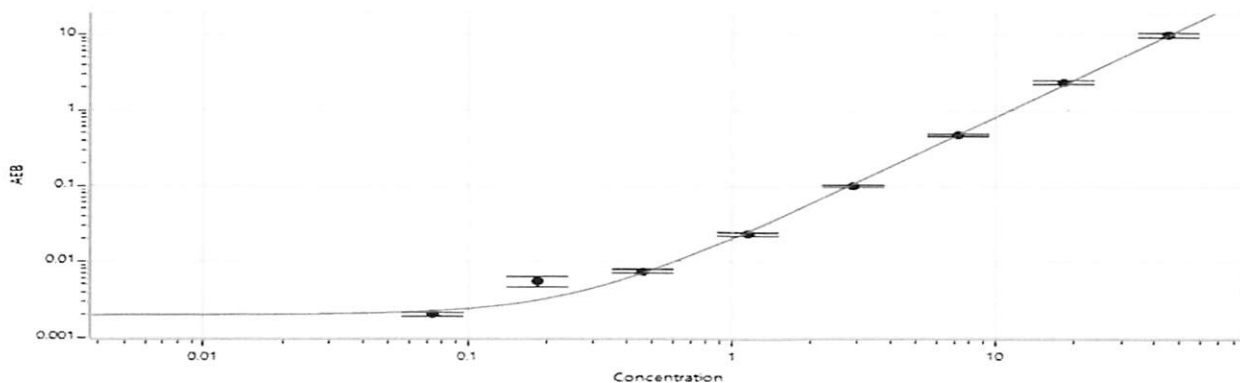
**Release Materials\***

**Abeta 40**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
<b>Control 1</b>	317329A	5.35	4.04 - 7.50
<b>Control 2</b>	317330A	89.8	61.6 - 114

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	972
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	972
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	972

Calibration Curve



**Calibrator Levels (pg/mL)**

A	0.000
B	0.074
C	0.185
D	0.463
E	1.16
F	2.89
G	7.21
H	18.1
I	45.2

**Release Materials\***

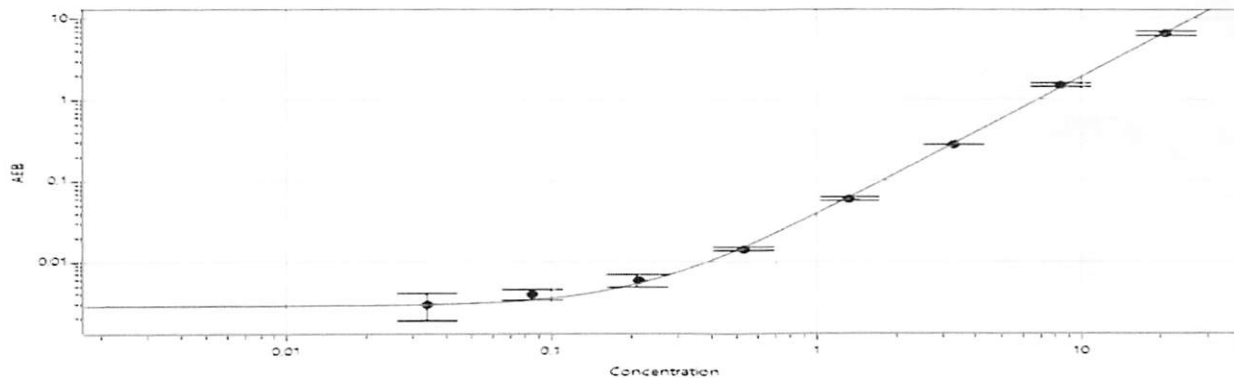
**Abeta 42**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
<b>Control 1</b>	317329A	2.37	1.83 - 3.40
<b>Control 2</b>	317330A	37.2	26.2 - 48.6

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

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	972
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	972
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	972

Calibration Curve



**Calibrator Levels (pg/mL)**

A	0.000
B	0.034
C	0.084
D	0.212
E	0.530
F	1.32
G	3.31
H	8.30
I	20.7

**Release Materials\***

**GFAP®**

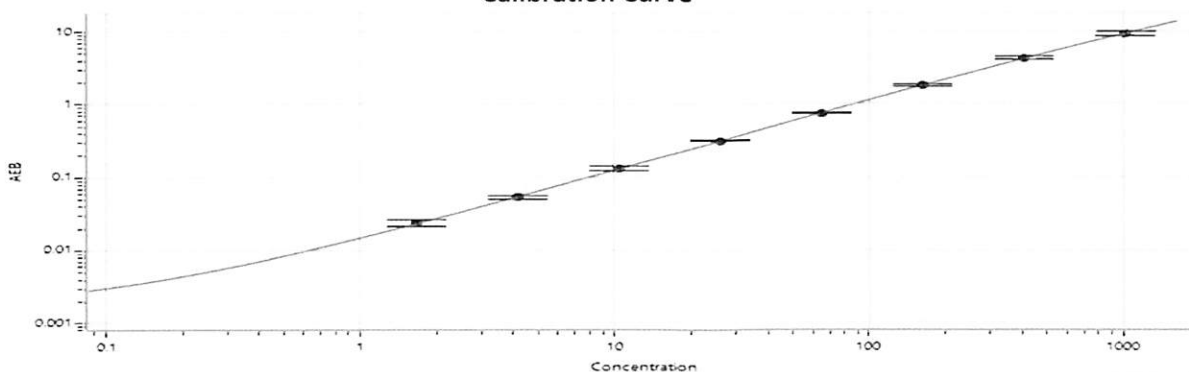
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
<b>Control 1</b>	317329A	122	88.6 - 190
<b>Control 2</b>	317330A	1778	1288 - 2392

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	972
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	972
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	972

**Calibrator Levels (pg/mL)**

<b>A</b>	0.000
<b>B</b>	1.67
<b>C</b>	4.18
<b>D</b>	10.4
<b>E</b>	26.1
<b>F</b>	65.3
<b>G</b>	163
<b>H</b>	408
<b>I</b>	1020

**Calibration Curve**



**Release Materials\***

**NF-Light®**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
<b>Control 1</b>	317329A	97.6	64.8 - 120
<b>Control 2</b>	317330A	1279	867 - 1610

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

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	972
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	972
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	972

**Calibrator Levels (pg/mL)**

<b>A</b>	0.000
<b>B</b>	1.10
<b>C</b>	2.75
<b>D</b>	6.88
<b>E</b>	17.2
<b>F</b>	43.0
<b>G</b>	107
<b>H</b>	269
<b>I</b>	672

**Calibration Curve**

