

Product Number:	103670
Lot Number:	503990
Expiration:	5-Oct-2024
Platform(s):	HD-X/HD-1

Component	Part Number	Lot Number
Bead Reagent	103653	327812
Detector Reagent	103655	327904
SBG Reagent	103657	327814
Plasma Sample Diluent	103659	327906
CSF Sample Diluent	103727	328211
Calibrators	103667	324402
Control 1	103668	324403
Control 2	103669	324403
RGP Reagent	103159	N/A <sup>1</sup>

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



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

Review/Approval

Cody Carlson QC Manager  
Name Title

*[Handwritten Signature]* 27 Dec 2023  
Signature/Date

**Release Materials\***

**Abeta 40**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	324403	18.1	15.3 - 23.0
Control 2	324403	85.1	80.9 - 133
Panel 1	226304	26.1	20.8 - 31.2
Panel 2	332403	55.1	44.4 - 68.3
Panel 3	333807	129	99.7 - 150

**Abeta 42**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	324403	8.22	6.72 - 10.1
Control 2	324403	36.4	30.8 - 46.1
Panel 1	226304	2.29	1.65 - 2.97
Panel 2	332403	10.6	7.78 - 11.7
Panel 3	333807	50.6	41.1 - 61.6

**GFAP®**

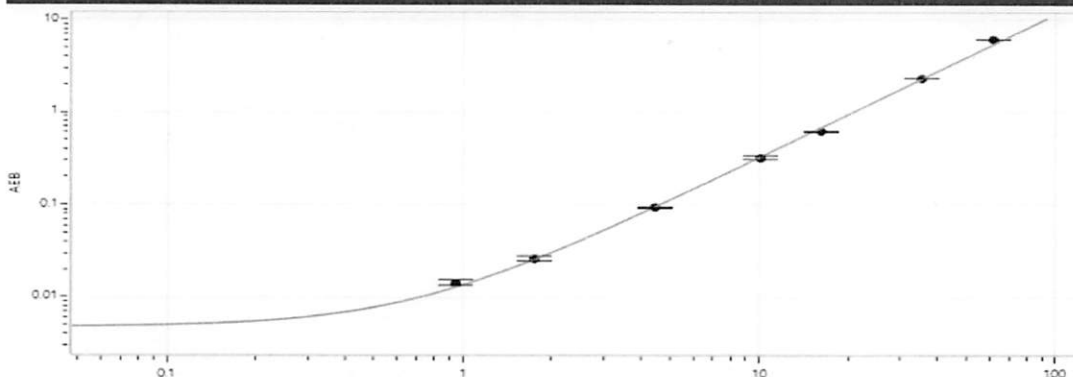
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	324403	171	149 - 223
Control 2	324403	3501	2955 - 4432
Panel 1	226304	39.8	28.1 - 58.0
Panel 2	332403	204	160 - 273
Panel 3	333807	911	728 - 1092

**NF-Light®**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	324403	22.2	19.0 - 28.5
Control 2	324403	495	419 - 629
Panel 1	226304	2.36	1.38 - 4.84
Panel 2	332403	64.4	59.5 - 89.2
Panel 3	333807	465	354 - 670

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

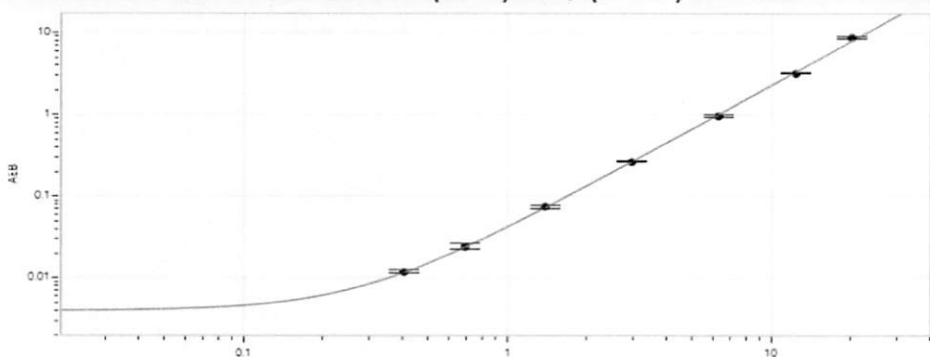
**Abeta 40 Calibrator Curve**



**Calibrator Levels  
(pg/mL)**

A	0.000
B	0.947
C	1.75
D	4.46
E	10.1
F	16.2
G	35.2
H	61.4

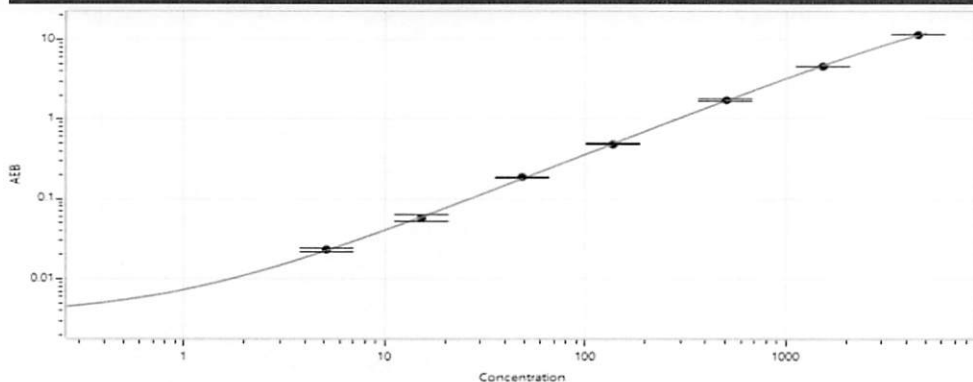
**Abeta 42 Calibrator Curve**



**Calibrator Levels  
(pg/mL)**

A	0.000
B	0.405
C	0.692
D	1.39
E	2.97
F	6.31
G	12.4
H	20.1

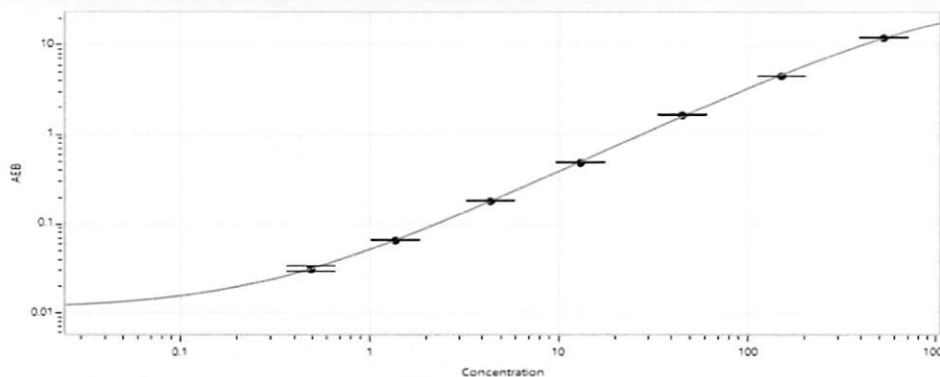
**GFAP® Calibrator Curve**



**Calibrator Levels  
(pg/mL)**

A	0.000
B	5.13
C	15.2
D	48.3
E	138
F	501
G	1529
H	4577

**NF-Light® Calibrator Curve**



**Calibrator Levels  
(pg/mL)**

A	0.000
B	0.489
C	1.38
D	4.38
E	13.0
F	45.5
G	152
H	525