

Product Number:	103670
Lot Number:	504173
Expiration:	5-Oct-2024
Platform(s):	HD-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	103653	327812	N/A
Detector Reagent	103655	327904	N/A
SBG Reagent	103657	327814	327814A
Plasma Sample Diluent	103659	327906	N/A
CSF Sample Diluent	103727	328211	N/A
Calibrators	103667	329705	N/A
Control 1	103668	329706	N/A
Control 2	103669	329706	N/A
RGP Reagent	103159	N/A ¹	N/A ¹

¹ Reagents are not Kit Lot Specific

Review/Approval

Cody Carlson
Name

QC Manager
Title

[Signature] 14 MAR 2024
Signature/Date

Release Materials*

Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	329706	16.7	14.8 - 22.2
Control 2	329706	102	100 - 150
Panel 1	226304	25.3	20.8 - 31.2
Panel 2	332403	60.9	45.3 - 67.9
Panel 3	401902	205	155 - 232

Abeta 42

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	329706	7.31	6.08 - 9.30
Control 2	329706	36.8	30.6 - 45.9
Panel 1	226304	2.31	1.65 - 2.97
Panel 2	332403	9.82	7.79 - 11.7
Panel 3	401902	73.2	55.2 - 82.8

GFAP®

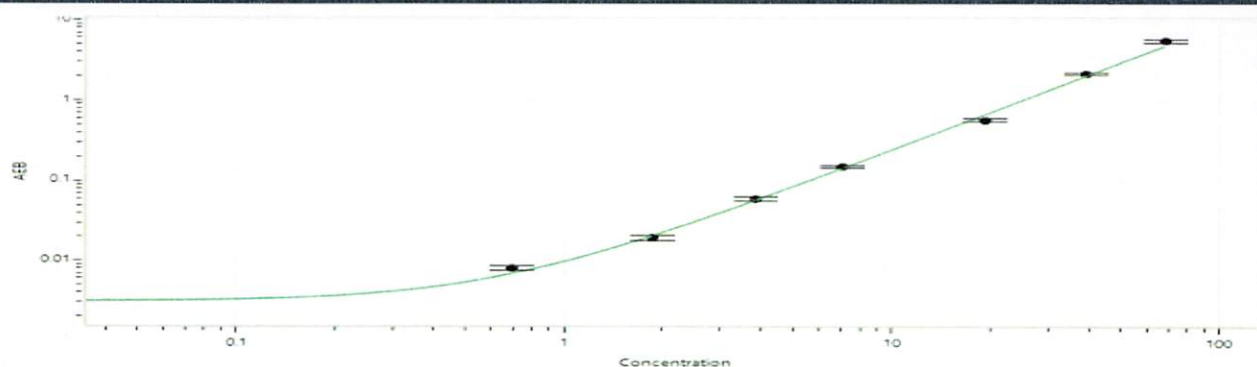
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	329706	183	166 - 248
Control 2	329706	4178	3507 - 5260
Panel 1	226304	42.4	28.1 - 58.0
Panel 2	332403	218	155 - 271
Panel 3	401902	4072	2497 - 4568

NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	329706	23.1	19.7 - 29.5
Control 2	329706	554	474 - 711
Panel 1	226304	2.87	1.38 - 4.84
Panel 2	332403	85.1	59.3 - 89.0
Panel 3	401902	705	367 - 769

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

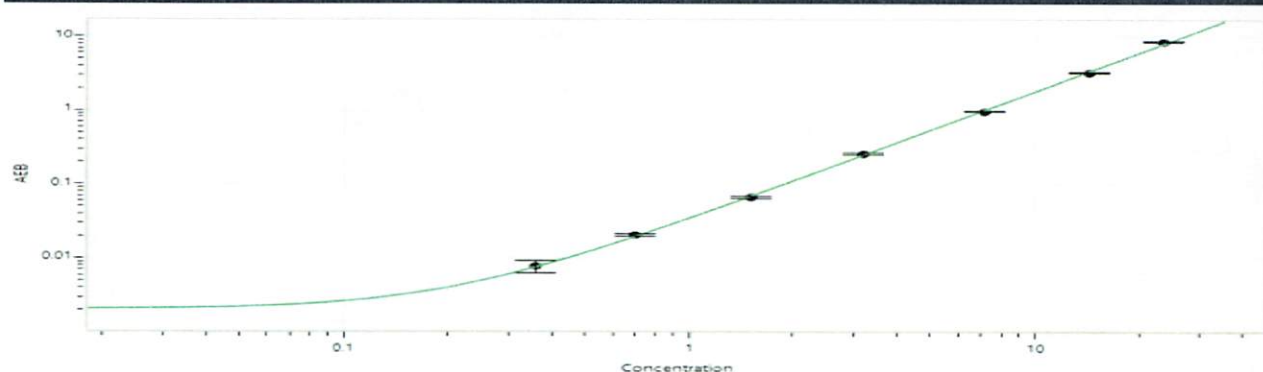
Abeta 40 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.693
C	1.87
D	3.83
E	7.07
F	19.2
G	39.4
H	68.3

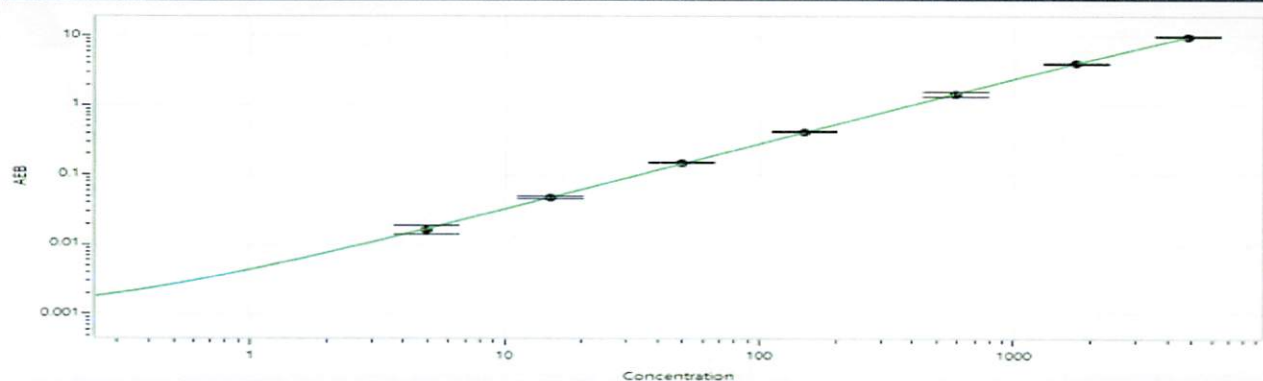
Abeta 42 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.360
C	0.700
D	1.50
E	3.20
F	7.11
G	14.3
H	23.5

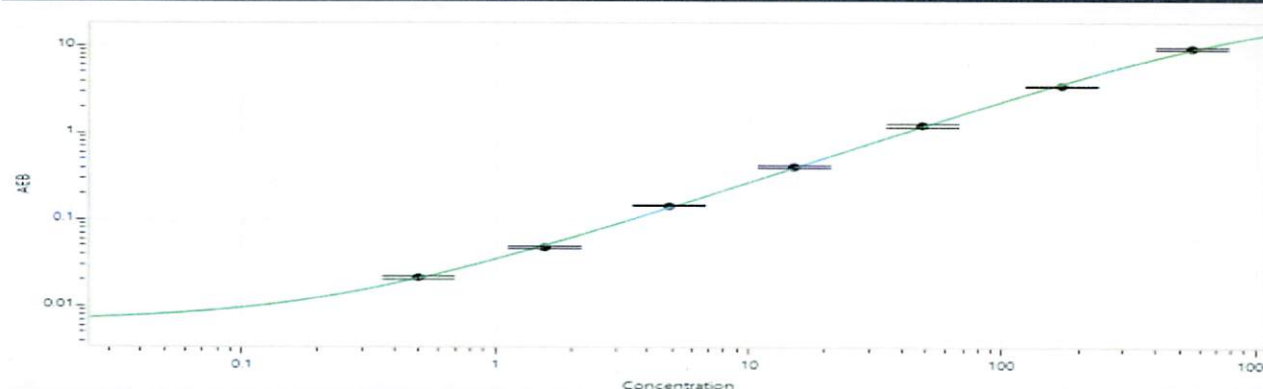
GFAP® Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	4.88
C	15.2
D	49.8
E	151
F	591
G	1757
H	4803

NF-Light® Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.498
C	1.56
D	4.82
E	15.0
F	48.3
G	170
H	558