

Product Number:	103670
Lot Number:	504166
Expiration:	5-Apr-2025
Platform(s):	HD-X



*Data below
represents results
generated on the*

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	324402	N/A
Control 1	103668	324403	N/A
Control 2	103669	324403	N/A
RGP Reagent	103159	N/A ¹	N/A ¹

¹ Reagents are not Kit Lot Specific

Review/Approval

Cody Carlson

QC Manager

Name

Title

 14 MAR 2024
Signature/Date

Release Materials*

Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	324403	17.1	15.3 - 23.0
Control 2	324403	94.7	80.9 - 133
Panel 1	226304	25.7	20.8 - 31.2
Panel 2	332403	59.8	45.3 - 67.9
Panel 3	401902	200	152 - 228

Abeta 42

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	324403	7.38	6.72 - 10.1
Control 2	324403	36.8	30.8 - 46.1
Panel 1	226304	2.20	1.65 - 2.97
Panel 2	332403	9.58	7.79 - 11.7
Panel 3	401902	68.9	54.4 - 81.6

GFAP®

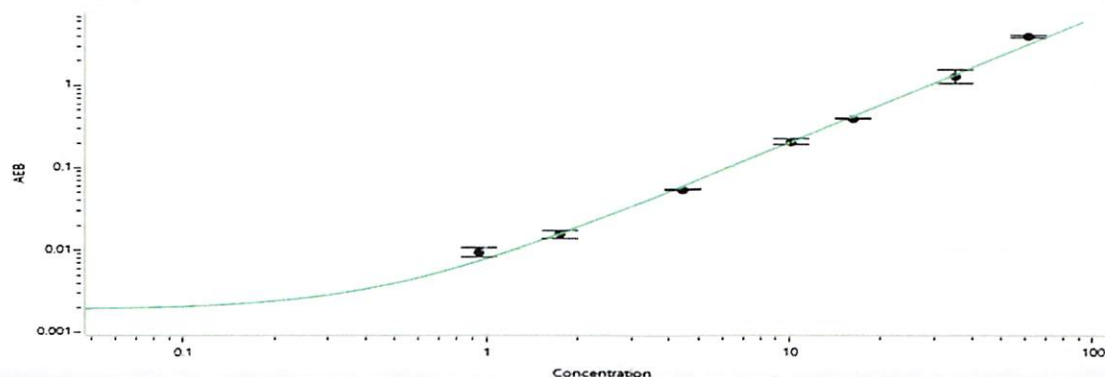
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	324403	159	149 - 223
Control 2	324403	3493	2955 - 4432
Panel 1	226304	39.0	28.1 - 58.0
Panel 2	332403	189	155 - 271
Panel 3	401902	3518	2529 - 4266

NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	324403	20.8	19.0 - 28.5
Control 2	324403	502	419 - 629
Panel 1	226304	2.61	1.38 - 4.84
Panel 2	332403	71.7	59.3 - 89.0
Panel 3	401902	575	423 - 645

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

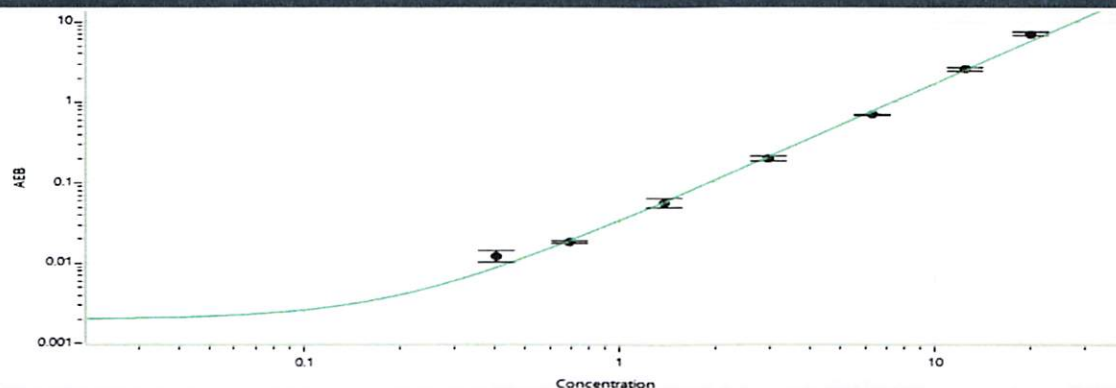
Abeta 40 Calibrator Curve



**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

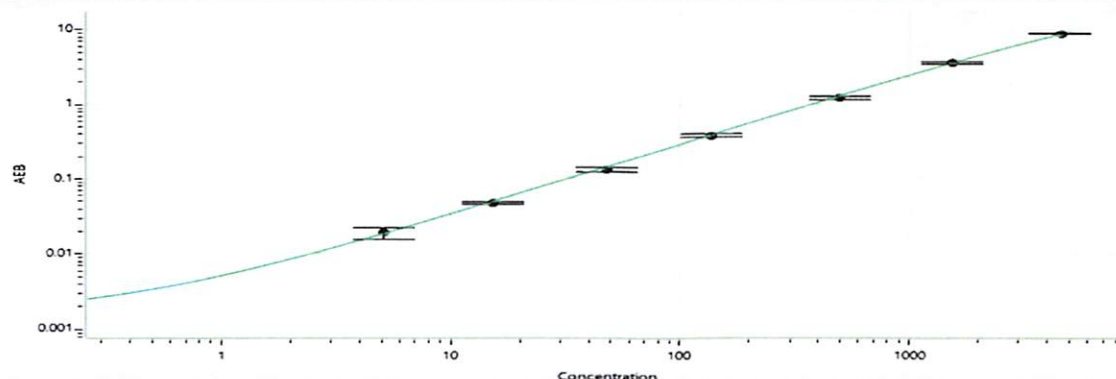
Abeta 42 Calibrator Curve



**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

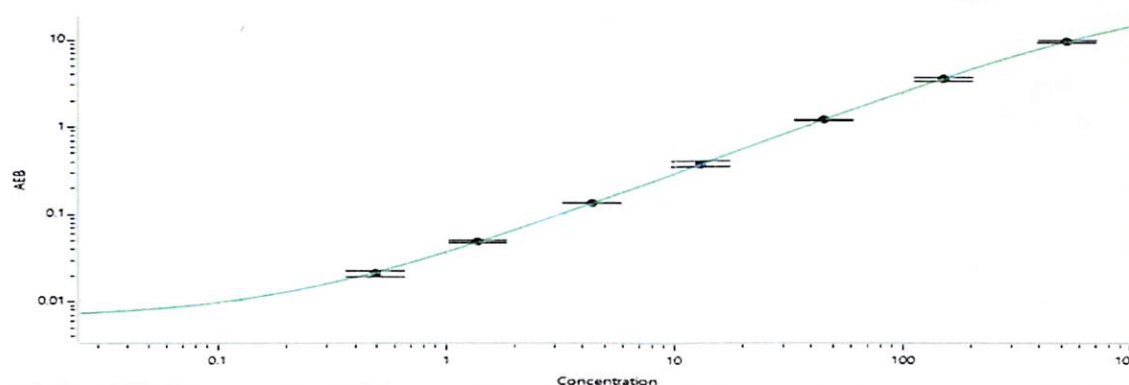
GFAP® Calibrator Curve



**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