

Product Number:	103670
Lot Number:	504197
Expiration:	8-Jun-2025
Platform(s):	HD-X

Data below represents results generated on the



Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	103653	334203	N/A
Detector Reagent	103655	334204	N/A
SBG Reagent	103657	334504	334504A
Plasma Sample Diluent	103659	334506	N/A
CSF Sample Diluent	103727	334505	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

QC Manager

Cody Carlson 10 APR 2024

Name

Title

Signature/Date

Release Materials*

Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	329706	17.1	14.6 - 21.9
Control 2	329706	116	98.7 - 148
Panel 1	226304	25.2	20.8 - 31.2
Panel 2	332403	55.5	45.1 - 67.7
Panel 3	401902	185	154 - 230

Abeta 42

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	329706	7.37	6.10 - 9.15
Control 2	329706	36.6	30.3 - 45.5
Panel 1	226304	1.92	1.65 - 2.97
Panel 2	332403	9.70	7.79 - 11.7
Panel 3	401902	68.3	55.1 - 82.7

GFAP®

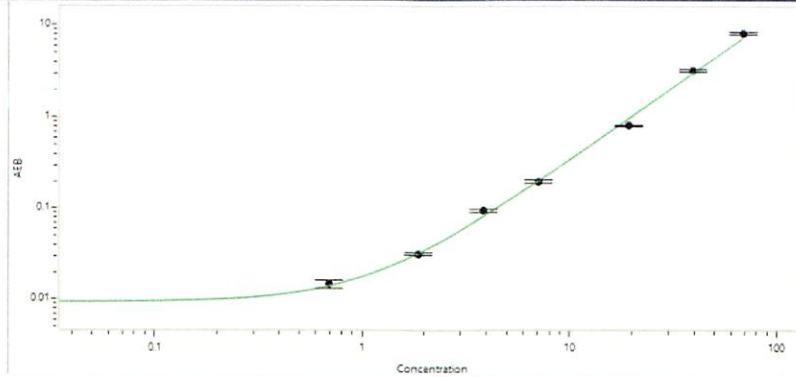
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	329706	181	161 - 242
Control 2	329706	3964	3440 - 5159
Panel 1	226304	37.8	28.1 - 58.0
Panel 2	332403	208	158 - 267
Panel 3	401902	3514	2584 - 4475

NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	329706	24.2	19.6 - 29.4
Control 2	329706	519	455 - 701
Panel 1	226304	2.63	1.38 - 4.84
Panel 2	332403	74.2	59.3 - 89.0
Panel 3	401902	576	385 - 753

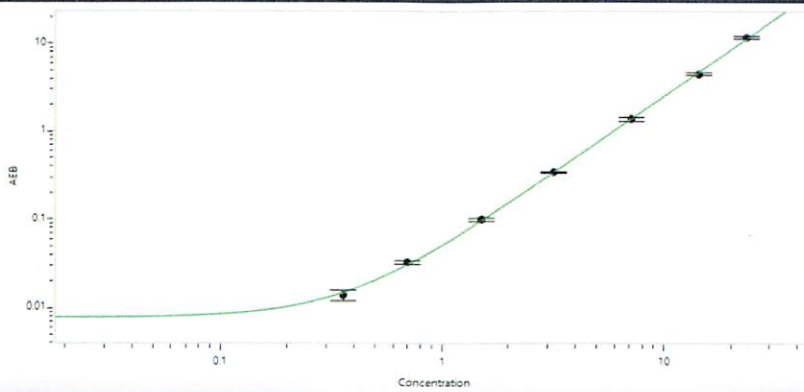
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Abeta 40 Calibrator Curve



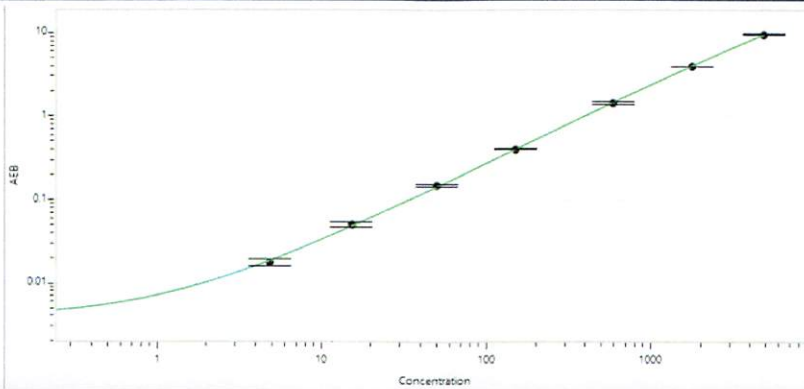
Calibrator Levels (pg/mL)	
A	0.000
B	0.693
C	1.87
D	3.83
E	7.07
F	19.2
G	39.4
H	68.3

Abeta 42 Calibrator Curve



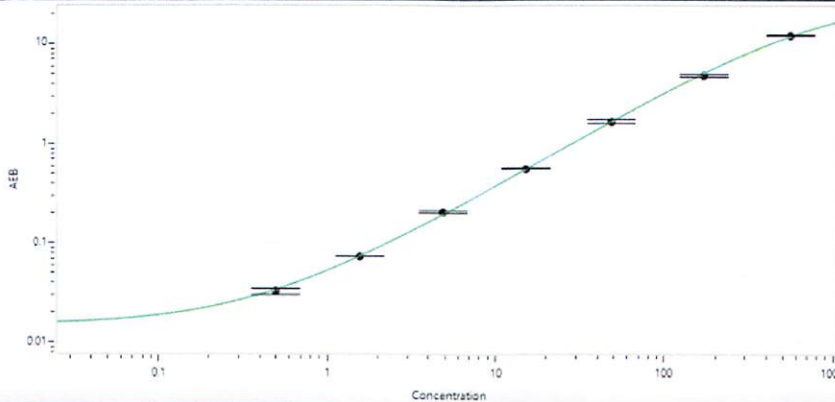
Calibrator Levels (pg/mL)	
A	0.000
B	0.360
C	0.700
D	1.50
E	3.20
F	7.11
G	14.3
H	23.5

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