

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

¹ Reagents are not Kit Lot Specific

Review/Approval

Dipesh Patel
Name

Sz. Doc. content - QA
Title

Signature/Date

Release Materials*
Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	329706	17.2	14.6 - 21.9
Control 2	329706	127	89.8 - 149

Abeta 42

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	329706	7.79	6.12 - 9.18
Control 2	329706	38.1	30.3 - 45.4

GFAP®

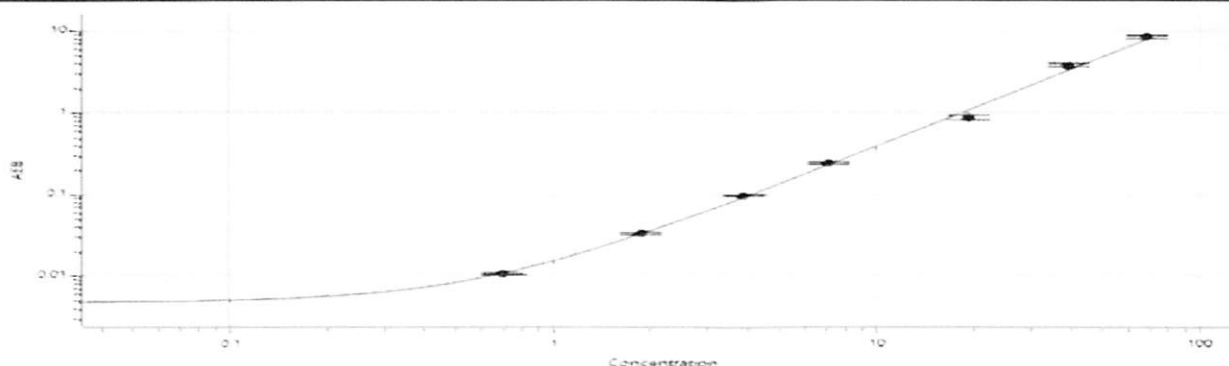
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	329706	202	161 - 242
Control 2	329706	4662	3450 - 5174

NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	329706	24.8	19.6 - 29.5
Control 2	329706	589	459 - 691

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

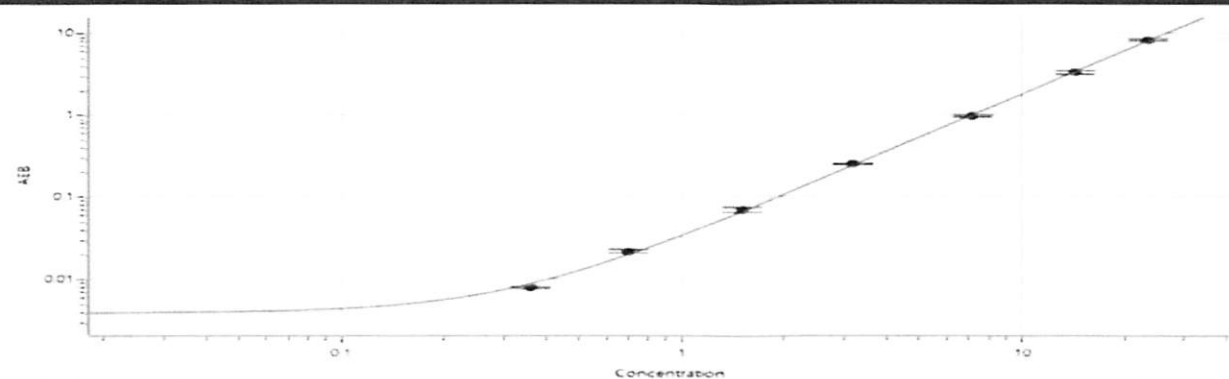
Abeta 40 Calibrator Curve



**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

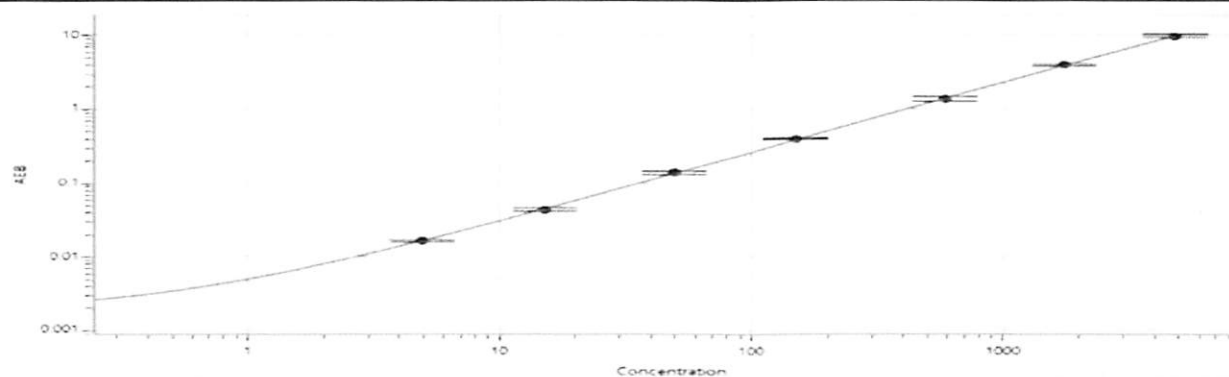
Abeta 42 Calibrator Curve



**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

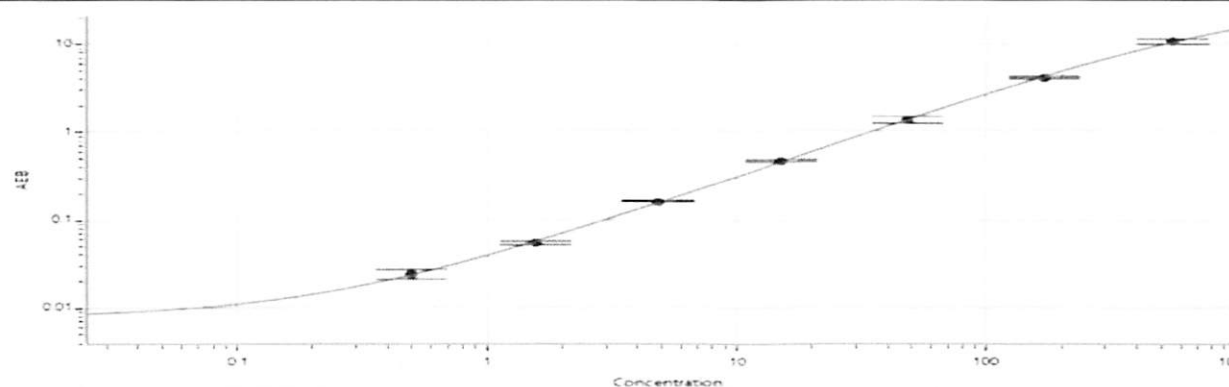
GFAP® Calibrator Curve



**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