

Product Number:	103670
Lot Number:	504582
Expiration:	26-Sep-2026
Platform(s):	HD-X

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	103653	511101	N/A
Detector Reagent	103655	511117	N/A
SBG Reagent	103657	511118	N/A
Plasma Sample Diluent	103659	511119	N/A
CSF Sample Diluent	103727	511116	511116A
Calibrators	103667	511106	N/A
Control 1	103668	511107	N/A
Control 2	103669	511107	N/A
RGP Reagent	103159	N/A ¹	N/A ¹

¹ Reagents are not Kit Lot Specific



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

Review/Approval

Dipesh Patel
Name

Sr. Doc. Control
Title

[Signature] | 12 MAY 2025
Signature/Date

Release Materials*
Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	511107	21.6	17.4 - 26.1
Control 2	511107	132	114 - 171

Abeta 42

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	511107	8.64	7.02 - 10.5
Control 2	511107	44.0	37.3 - 55.9

GFAP®

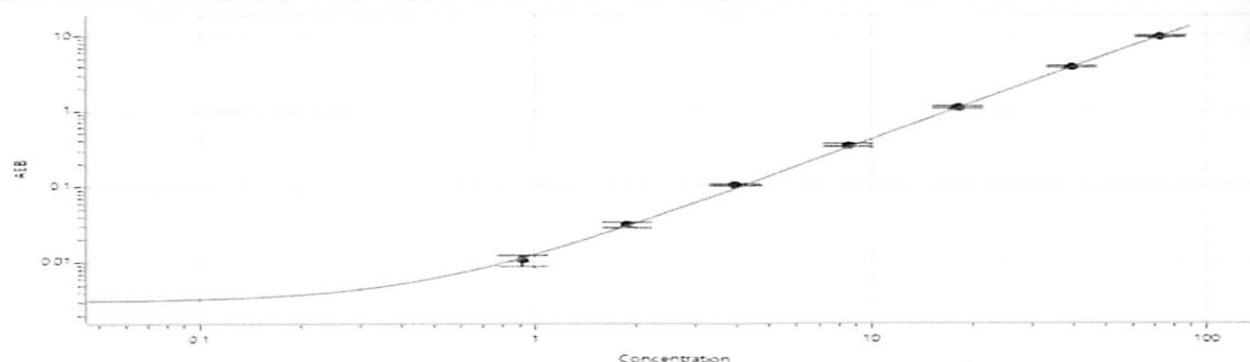
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	511107	214	179 - 268
Control 2	511107	3682	3383 - 5532

NF-Light®

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	511107	24.6	20.2 - 30.4
Control 2	511107	477	388 - 669

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

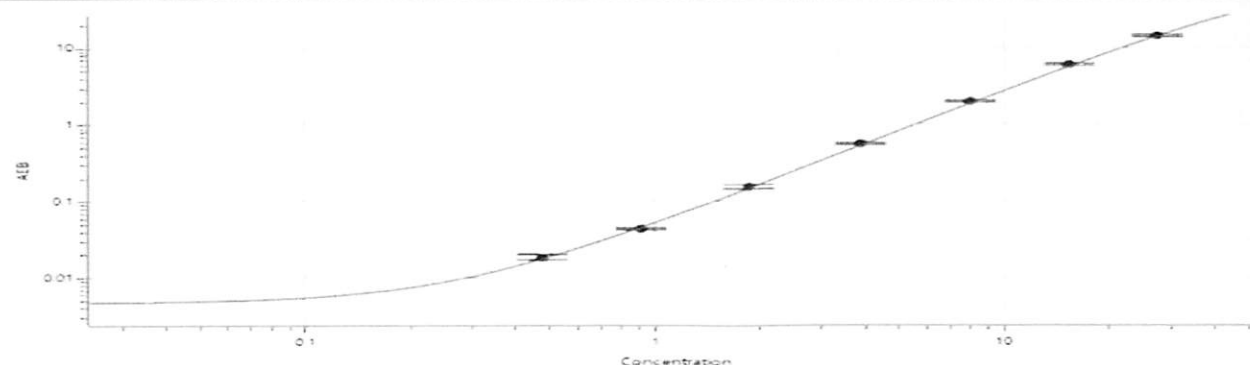
Abeta 40 Calibrator Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.914
C	1.88
D	3.93
E	8.57
F	18.1
G	39.5
H	72.9

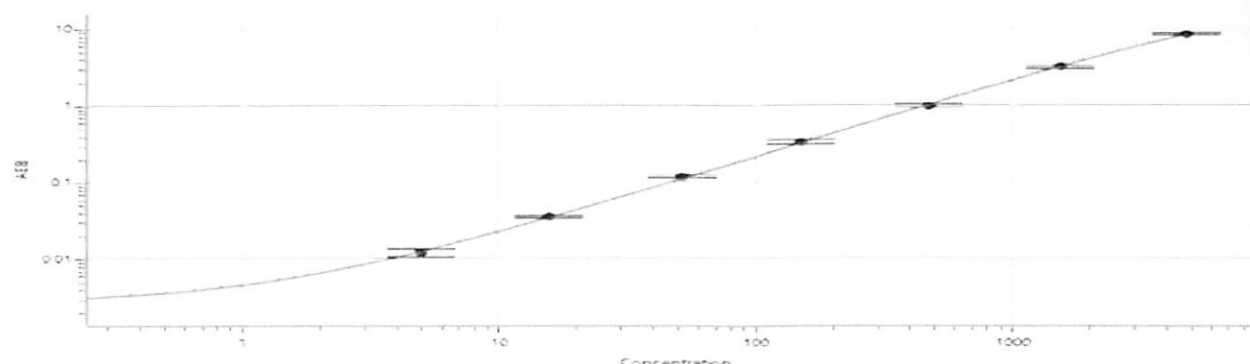
Abeta 42 Calibrator Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.477
C	0.908
D	1.85
E	3.86
F	8.02
G	15.3
H	27.4

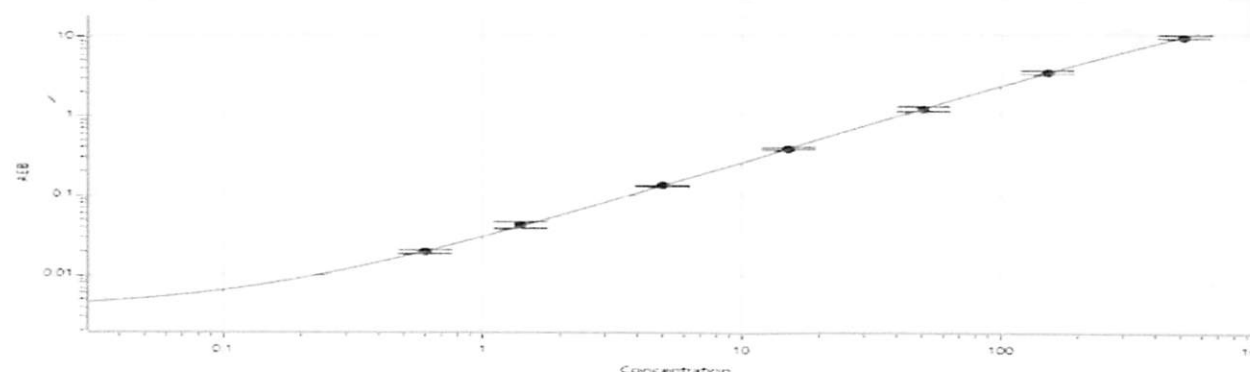
GFAP® Calibrator Curve



Calibrator Levels (pg/mL)

A	0.000
B	4.94
C	15.7
D	52.1
E	151
F	478
G	1539
H	4767

NF-Light® Calibrator Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.602
C	1.40
D	5.02
E	15.2
F	50.6
G	153
H	516