

Product Number:	104712
Lot Number:	504329
Expiration:	26-Jan-2026
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	104754	421402
Detector Reagent	104696	425413
SBG Reagent	104698	425705
Plasma Sample Diluent	104716	421215
CSF Sample Diluent	104717	421115
Calibrator Diluent	104699	420806
Control Diluent	104715	420807
Calibrator Concentrate	104708	403023
Control 1 Concentrate	104710	403024
Control 2 Concentrate	104711	403025
RGP 2.0 Reagent	104321	N/A [†]
Activation Buffer Reagent	104324	N/A [†]

[†] Reagents are not Kit Lot Specific

Review/Approval

Kyle Giroux
Name

Senior Document Control Specialist, QA
Title

[Signature] 29 Oct 2024
Signature/Date

Release Materials*

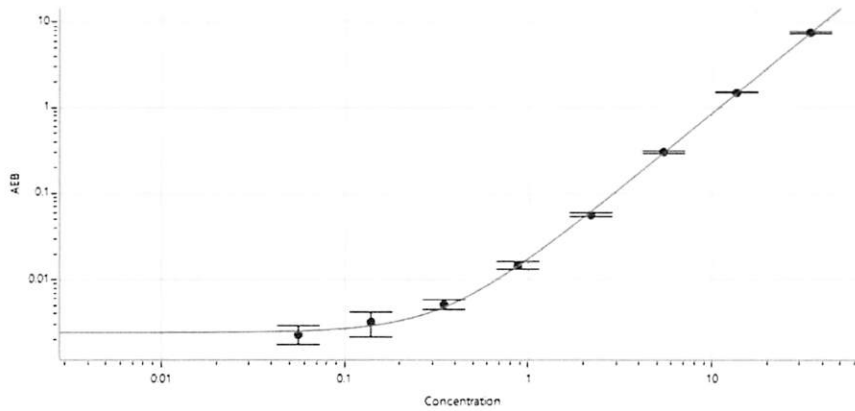
Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	403024	4.36	2.82 - 5.24
Control 2	403025	62.6	41.9 - 82.8

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

Calibrator Concentrate Reconstitution Volume (uL)	1144
Control 1 Concentrate Reconstitution Volume (uL)	1144
Control 2 Concentrate Reconstitution Volume (uL)	1144

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.056
C	0.140
D	0.349
E	0.872
F	2.19
G	5.45
H	13.6
I	34.1

Release Materials*

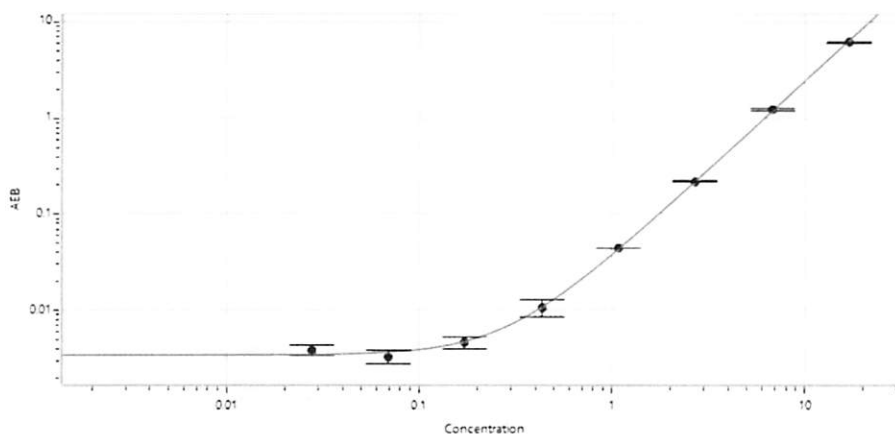
Abeta 42

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	403024	2.12	1.13 - 2.80
Control 2	403025	27.3	19.2 - 35.6

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

Calibrator Concentrate Reconstitution Volume (uL)	1144
Control 1 Concentrate Reconstitution Volume (uL)	1144
Control 2 Concentrate Reconstitution Volume (uL)	1144

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.028
C	0.069
D	0.173
E	0.432
F	1.08
G	2.70
H	6.75
I	16.8