

Component	Part Number	Lot Number
Bead Reagent	104754	510801
Detector Reagent	104696	425413
SBG Reagent	104698	425705
Plasma Sample Diluent	104716	422525
CSF Sample Diluent	104717	428510
Calibrator Diluent	104699	426819
Control Diluent	104715	426820
Calibrator Concentrate	104708	331021
Control 1 Concentrate	104710	403024
Control 2 Concentrate	104711	403025
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

Review/Approval

Name: Dipesh Patel Title: Sr. Doc. control -QA Signature/Date: [Signature] / 23 May 2025

Release Materials*

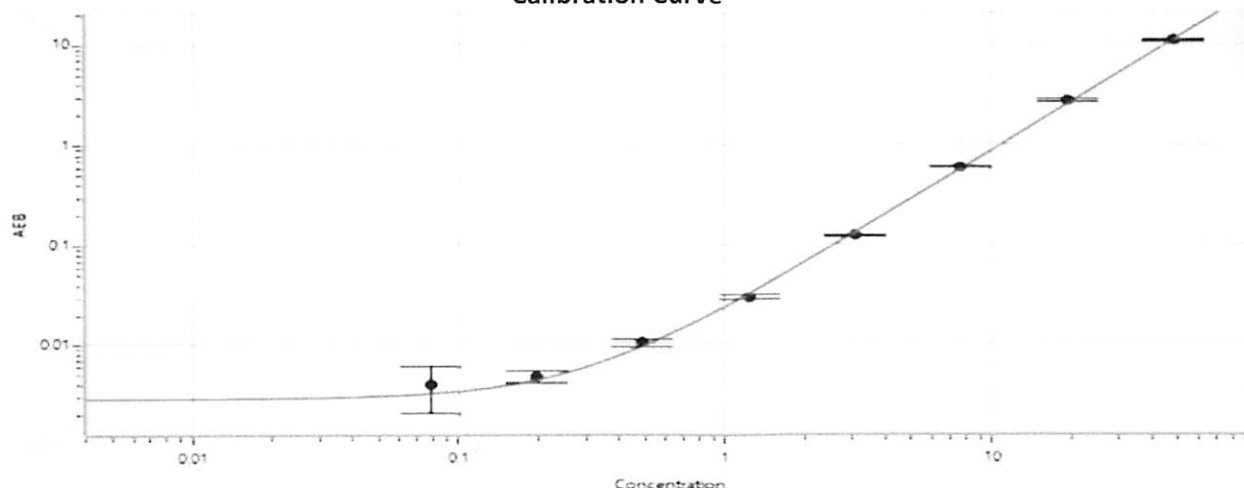
Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	403024	5.56	3.38 - 6.28
Control 2	403025	83.5	57.9 - 107

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

Calibrator Concentrate Reconstitution Volume (uL)	972
Control 1 Concentrate Reconstitution Volume (uL)	972
Control 2 Concentrate Reconstitution Volume (uL)	972

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.079
C	0.196
D	0.491
E	1.23
F	3.07
G	7.65
H	19.2
I	47.9

Release Materials*

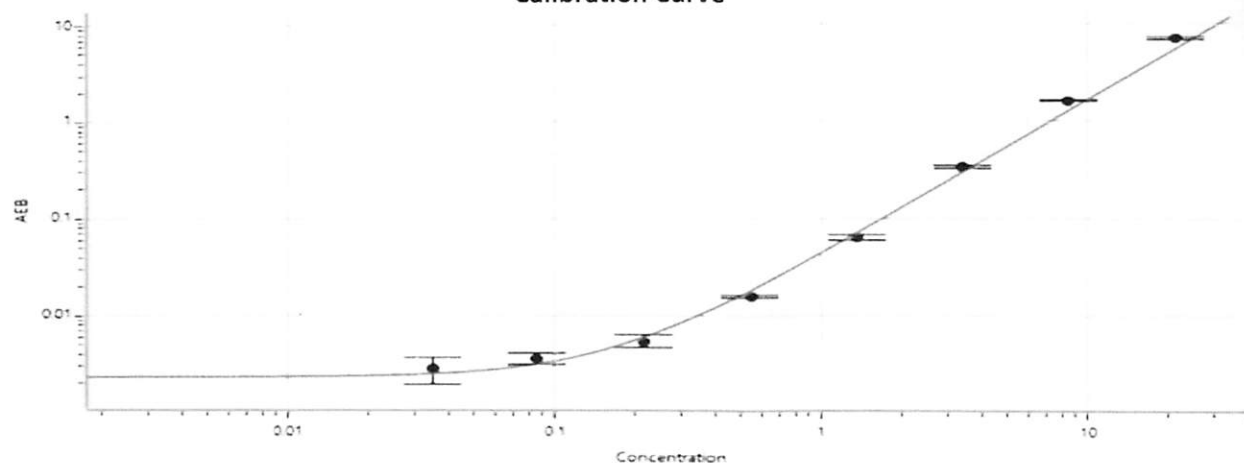
Abeta 42

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	403024	2.26	1.91 - 3.59
Control 2	403025	41.8	26.4 - 49.1

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

Calibrator Concentrate Reconstitution Volume (uL)	972
Control 1 Concentrate Reconstitution Volume (uL)	972
Control 2 Concentrate Reconstitution Volume (uL)	972

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.035
C	0.086
D	0.217
E	0.542
F	1.35
G	3.39
H	8.49
I	21.2