

Product Number:	104712
Lot Number:	504091
Expiration:	7-Jan-2025
Platform(s):	HD-X



Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	104754	334516	N/A
Detector Reagent	104696	334707	N/A
SBG Reagent	104698	334706	N/A
Plasma Sample Diluent	104716	329812	N/A
CSF Sample Diluent	104717	329625	N/A
Calibrator Diluent	104699	318813	318813A
Control Diluent	104715	318814	318814A
Calibrator Concentrate	104708	403015	N/A
Control 1 Concentrate	104710	403016	N/A
Control 2 Concentrate	104711	403017	N/A
RGP 2.0 Reagent	104321	N/A [†]	N/A [†]
Activation Buffer Reagent	104324	N/A [†]	N/A [†]

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

[†] Reagents are not Kit Lot Specific

Review/Approval

Cody Carlson
Name

QC Manager
Title

Cody Carlson 21 MAR 2024
Signature/Date

Release Materials*

Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	403016	3.95	2.87 - 5.33
Control 2	403017	64.6	44.7 - 83.0
Panel 1	327811	3.79	2.72 - 5.05
Panel 2	327811	39.8	30.0 - 55.7

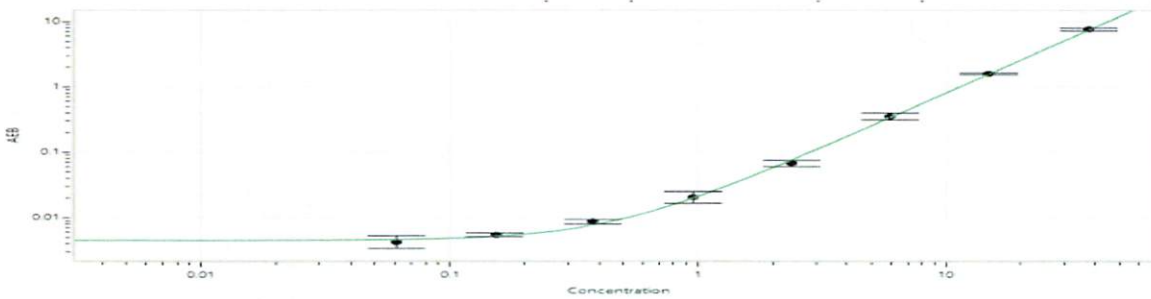
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

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

Calibrator Levels (pg/mL)

A	0.000
B	0.061
C	0.153
D	0.381
E	0.953
F	2.39
G	5.96
H	14.9
I	37.3

Calibration Curve



Release Materials*

Abeta 42

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	403016	2.20	1.66 - 3.09
Control 2	403017	35.7	25.5 - 47.4
Panel 1	327811	3.53	2.83 - 5.25
Panel 2	327811	33.7	24.4 - 45.4

* 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

Calibrator Levels (pg/mL)

A	0.000
B	0.039
C	0.097
D	0.243
E	0.607
F	1.51
G	3.79
H	9.48
I	23.6

Calibration Curve

