

Product Number:	101995
Lot Number:	504548
Expiration:	3-Sep-2026
Platform(s):	SR-X and HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	101998	506404
Detector Reagent	101996	506405
SBG Reagent	102000	506406
Sample Diluent	102002	506407
Reference Calibrators	103221	506907
Control 1	103222	506908
Control 2	103223	506908
SBG Diluent (SR-X) [†]	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

[†] Specific to use on the SR-X platform - refer to Neuro-3 Plex A Kit Instructions SR-X (KI-0083).

Review/Approval

Dhruvi Patel DOC Control II, QA SDP 27 Mar 2025

Name

Title

Signature/Date

Release Materials***Tau**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	506908	2.53	1.94 - 2.91
Control 2	506908	112	87.3 - 131

Abeta 42

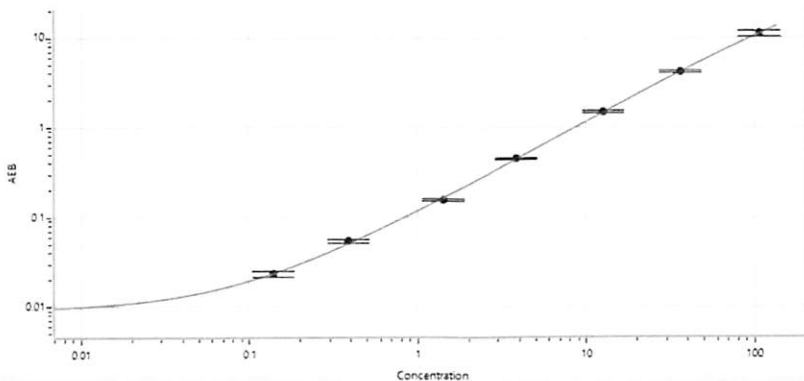
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	506908	3.70	2.78 - 4.40
Control 2	506908	94.4	75.9 - 114

Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	506908	9.67	5.25 - 11.0
Control 2	506908	158	121 - 181

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

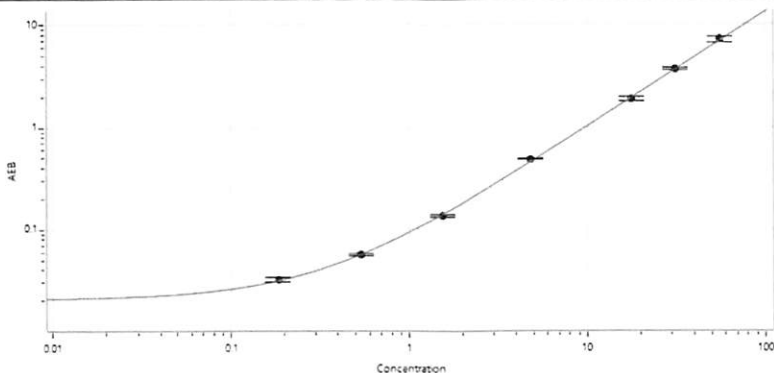
Tau Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.138
C	0.389
D	1.42
E	3.82
F	12.5
G	36.3
H	106

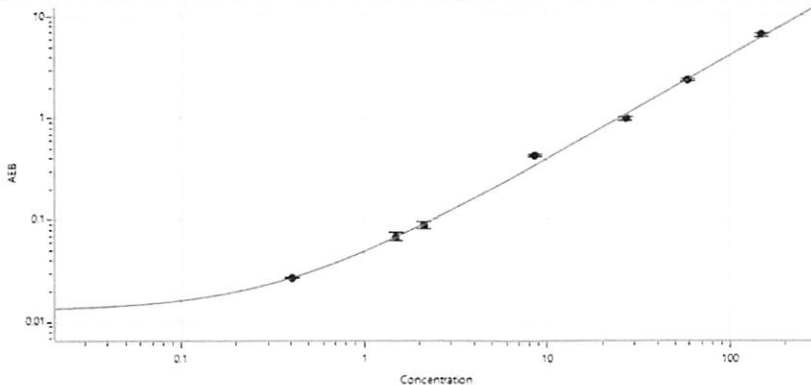
Abeta 42 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.184
C	0.535
D	1.53
E	4.83
F	17.7
G	31.4
H	55.8

Abeta 40 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.407
C	1.50
D	2.13
E	8.57
F	26.8
G	58.8
H	149