

Product Number:	101995
Lot Number:	504706
Expiration:	27-Feb-2026
Platform(s):	SR-X and HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	101998	523704
Detector Reagent	101996	523301
SBG Reagent	102000	523307
Sample Diluent	102002	523306
Reference Calibrators	103221	412804
Control 1	103222	405802
Control 2	103223	405802
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).



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

Review/Approval

Dhruvi Patel

Name

QA Specialist II, QA

Title

[Signature] *22SEP2025*

Signature/Date

Release Materials*

Tau

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	405802	2.26	1.87 - 2.81
Control 2	405802	112	92.6 - 139

Abeta 42

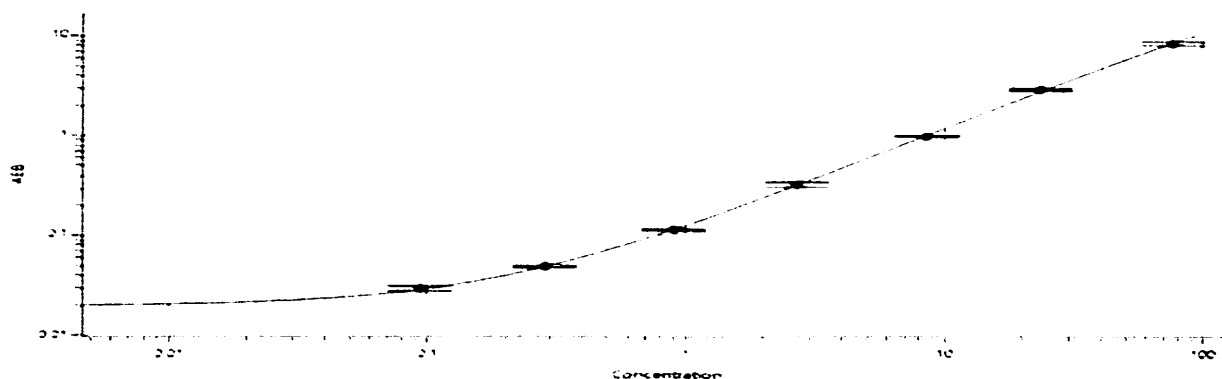
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	405802	3.13	2.63 - 3.95
Control 2	405802	82.6	70.8 - 106

Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	405802	9.31	6.70 - 10.1
Control 2	405802	163	140 - 210

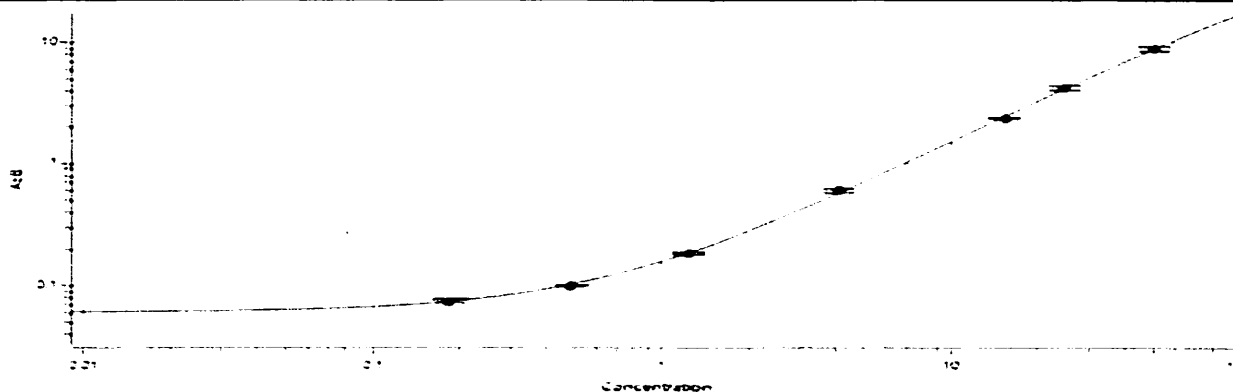
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Tau Calibrator Curve



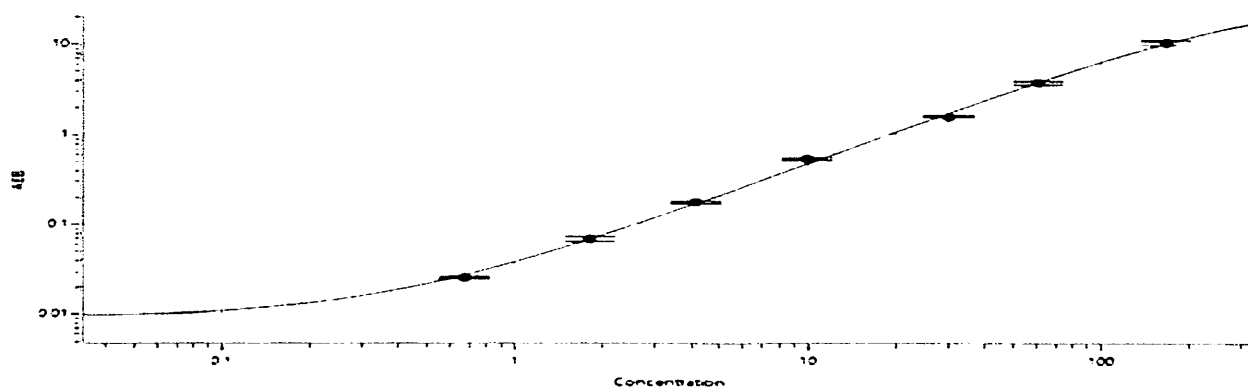
Calibrator Levels (pg/mL)	
A	0.000
B	0.093
C	0.285
D	0.902
E	2.69
F	8.50
G	23.5
H	76.0

Abeta 42 Calibrator Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.182
C	0.481
D	1.23
E	4.10
F	15.3
G	24.8
H	50.2

Abeta 40 Calibrator Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.671
C	1.80
D	4.11
E	9.85
F	30.1
G	61.0
H	167