

Product Number:	101995
Lot Number:	504214
Expiration:	2-Sep-2024
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	Control Lot Number
Bead Reagent	101998	306107	N/A
Detector Reagent	101996	306201	N/A
SBG Reagent	102000	306202	N/A
Sample Diluent	102002	306203	N/A
Calibrators	103221	317710	317710A
Control 1	103222	316407	316407A
Control 2	103223	316407	316407A
SBG Diluent (SRX)†	103158	N/A ¹	N/A ¹
RGP Reagent	103159	N/A ¹	N/A ¹

¹ Reagents are not Kit Lot Specific

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

Review/Approval

Cody Carlson

QC Manager

Name

Title

Cody Carlson 15 APR 2024
Signature/Date

Release Materials*

Tau

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	316407A	1.53	1.24 - 1.85
Control 2	316407A	63.2	53.2 - 79.7
Panel 1	124607	2.67	2.54 - 3.81
Panel 2	208403	126	86.2 - 177
Panel 3	208403	2.15	1.91 - 3.59

Abeta 42

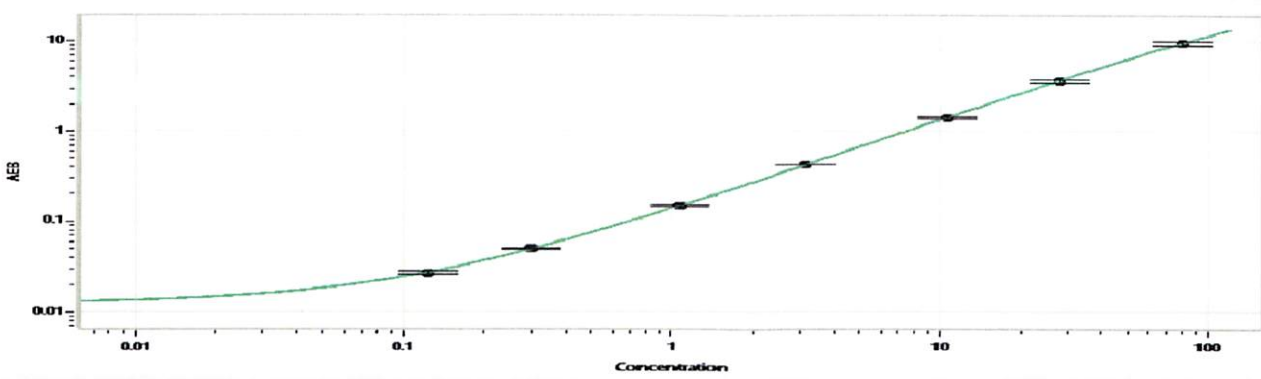
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	316407A	2.07	1.59 - 2.44
Control 2	316407A	80.5	60.8 - 91.1
Panel 1	124607	5.22	4.55 - 7.43
Panel 2	208403	7.18	6.55 - 10.3
Panel 3	208403	94.0	80.8 - 145

Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	316407A	5.42	3.87 - 6.08
Control 2	316407A	130	109 - 164
Panel 1	124607	123	114 - 171
Panel 2	208403	128	101 - 203
Panel 3	208403	78.7	63.8 - 118

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

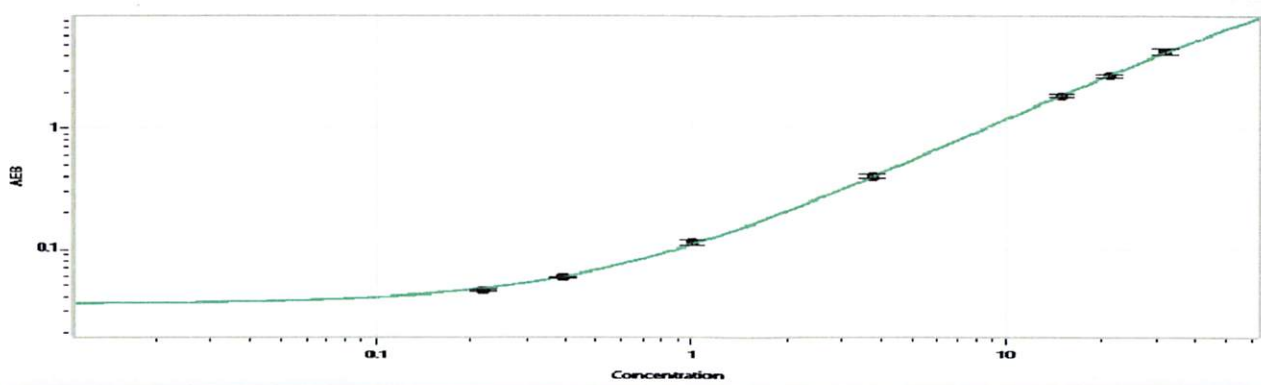
Tau Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.123
C	0.300
D	1.07
E	3.16
F	10.5
G	27.8
H	80.0

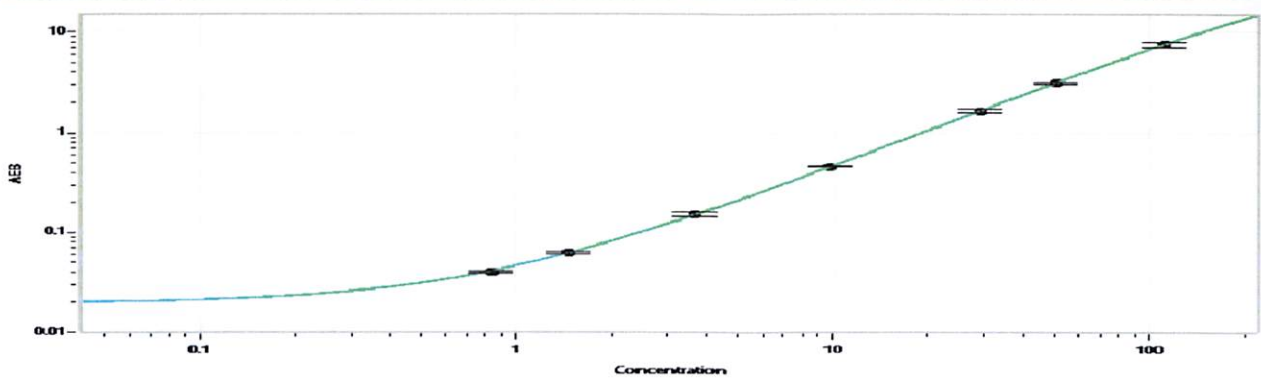
Abeta 42 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.219
C	0.390
D	1.01
E	3.75
F	14.9
G	21.2
H	31.9

Abeta 40 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.835
C	1.45
D	3.67
E	9.75
F	29.2
G	50.6
H	111