

Product Number:	101995
Lot Number:	504004
Expiration:	8-Nov-2024
Platform(s):	SR-X and HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	101998	333604
Detector Reagent	101996	333507
SBG Reagent	102000	333508
Sample Diluent	102002	333511
Calibrators	103221	333408
Control 1	103222	333409
Control 2	103223	333409
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

Cody Carlson
Name

QC Manager
Title

[Signature] 29 Dec 2023
Signature/Date

Release Materials*

Tau

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	333409	1.47	1.09 - 1.64
Control 2	333409	63.6	48.7 - 73.1
Panel 1	319110A	3.43	2.74 - 4.11
Panel 2	208403	138	110 - 167
Panel 3	208403	2.36	1.61 - 2.84

Abeta 42

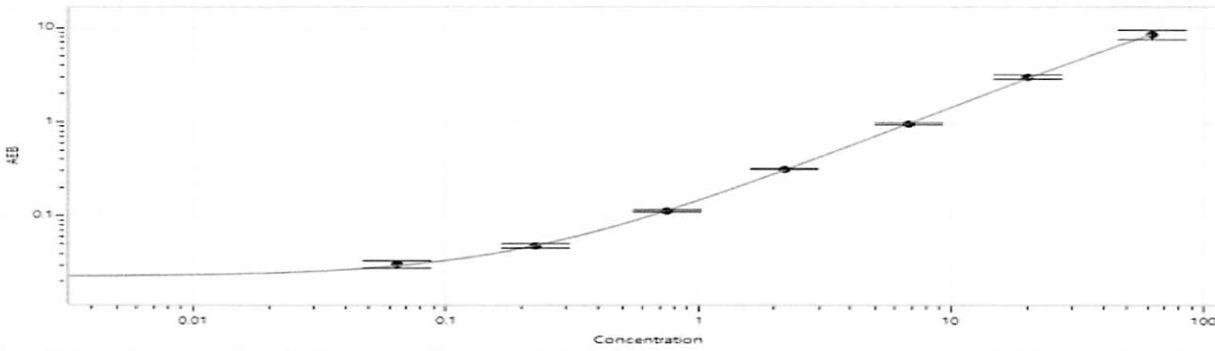
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	333409	4.14	2.42 - 4.68
Control 2	333409	97.6	68.5 - 112
Panel 1	319110A	2.92	2.30 - 3.45
Panel 2	208403	9.98	5.69 - 11.3
Panel 3	208403	132	80.4 - 155

Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	333409	11.9	7.38 - 13.1
Control 2	333409	172	142 - 213
Panel 1	319110A	78.9	63.4 - 95.0
Panel 2	208403	130	105 - 185
Panel 3	208403	79.0	60.6 - 113

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

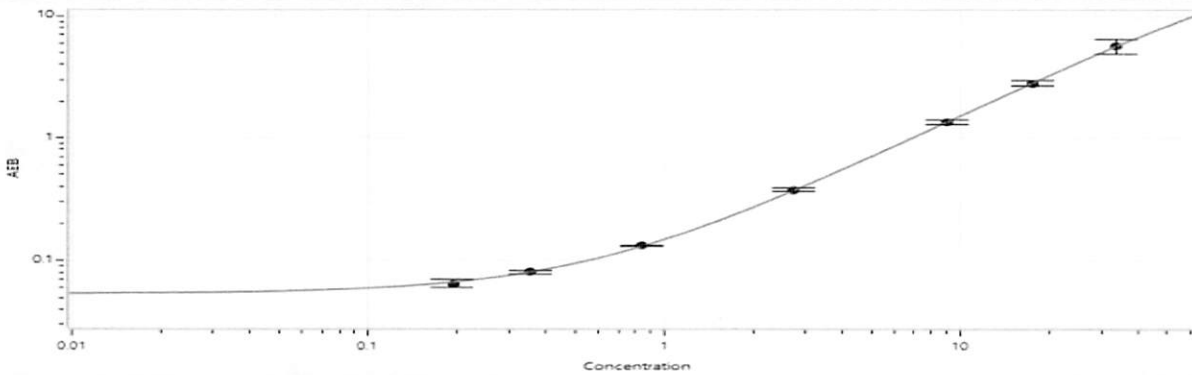
Tau Calibrator Curve



Calibrator Levels (pg/mL)

A	0
B	0.064
C	0.227
D	0.757
E	2.19
F	6.79
G	20.2
H	61.9

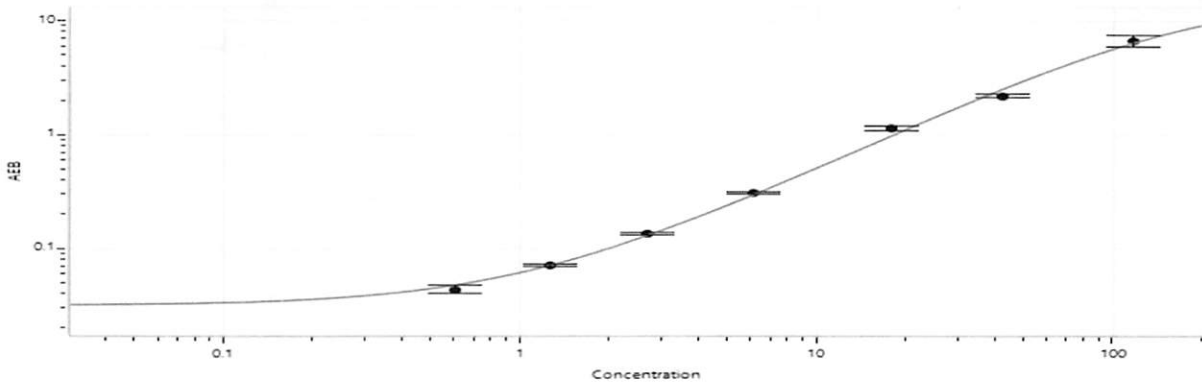
Abeta 42 Calibrator Curve



Calibrator Levels (pg/mL)

A	0
B	0.193
C	0.353
D	0.841
E	2.73
F	9.01
G	17.6
H	33.7

Abeta 40 Calibrator Curve



Calibrator Levels (pg/mL)

A	0
B	0.604
C	1.27
D	2.70
E	6.15
F	18.0
G	42.6
H	117