

Product Number:	101995
Lot Number:	503863
Expiration:	27-Jun-2024
Platform(s):	HD-1/HD-X & SR-X



Component	Part Number	Lot Number	Control Number
Bead Reagent	101998	317809	317809A
Detector Reagent	101996	318102	318102A
SBG Reagent	102000	318101	318101A
Sample Diluent	102002	317904	317904A
Calibrators	103221	317710	317710A
Control 1	103222	315103	315103A
Control 2	103223	315103	315103A
SBG Diluent (SRX)†	103158	N/A ¹	N/A ¹
RGP Reagent	103159	N/A ¹	N/A ¹

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

¹ Reagents are not Kit Lot Specific

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

Release Materials

Tau

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	315103	1.18	1.02 - 1.54
Control 2	315103	52.5	39.6 - 60.9
Panel 1	319110	3.03	2.74 - 4.11
Panel 2	208403	143	109 - 169
Panel 3	208403	1.94	1.57 - 2.82

Abeta 42

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	315103	2.36	1.67 - 3.23
Control 2	315103	57.4	39.5 - 76.7
Panel 1	319110	2.46	2.28 - 3.43
Panel 2	208403	8.38	5.86 - 10.6
Panel 3	208403	113	79.2 - 150

Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	315103	7.84	5.88 - 9.43
Control 2	315103	127	90.2 - 175
Panel 1	319110	73.7	63.4 - 95.2
Panel 2	208403	138	109 - 187
Panel 3	208403	76.1	62.0 - 115

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

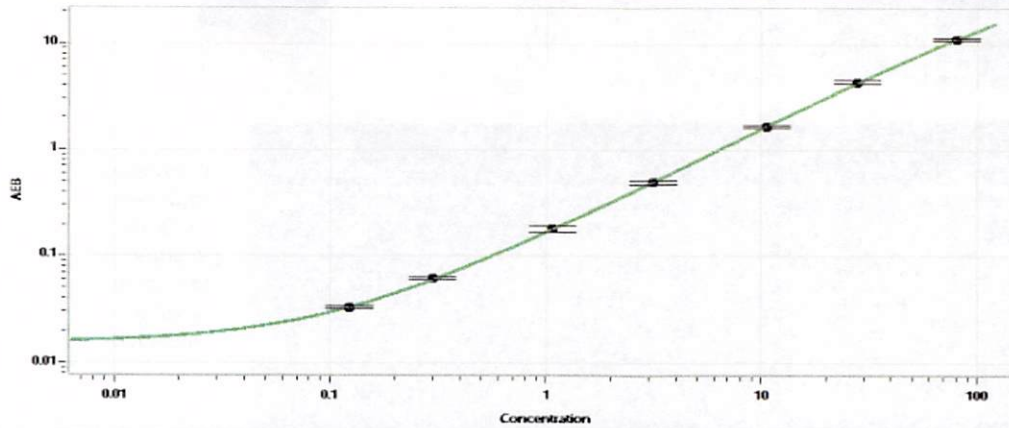
Review/Approval

Sonia Khillan
Name

QC
Title

Sonia Khillan 17AUG2023
Signature/Date

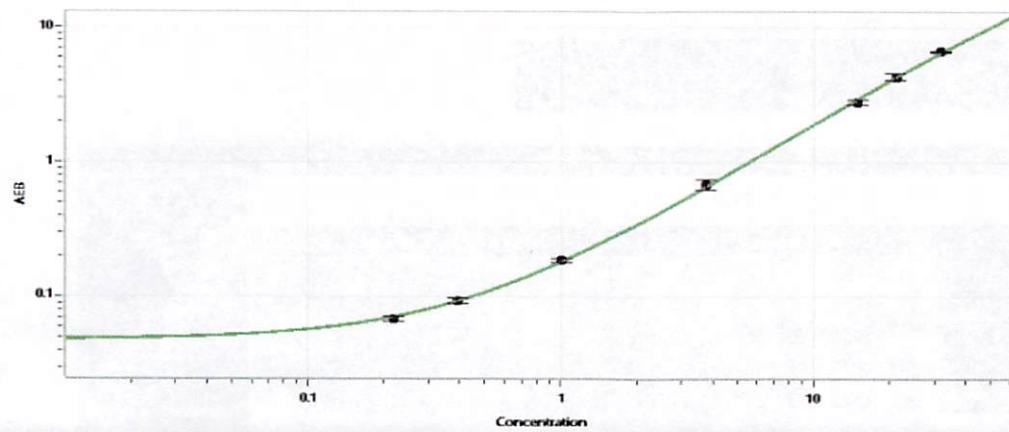
Tau Calibration Curve



**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

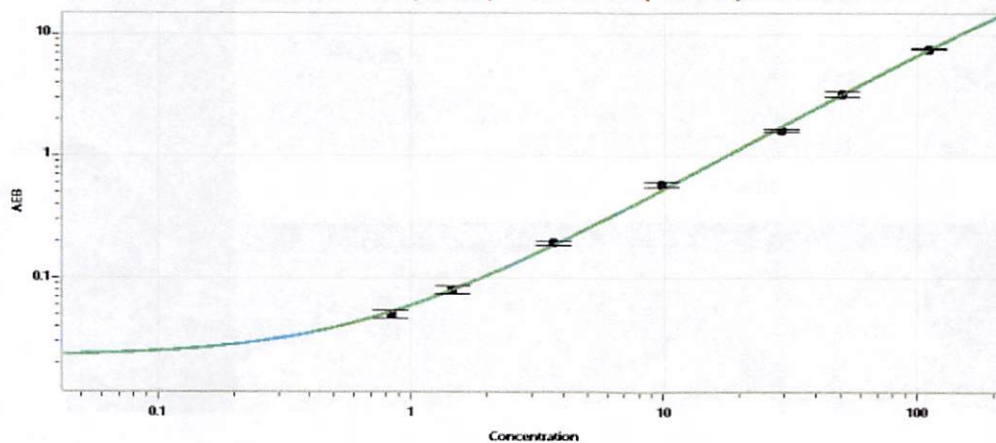
Abeta 42 Calibration Curve



**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 Calibration 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