

Product Number:	101995
Lot Number:	503659
Expiration:	8-Sep-2023
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101998	230707
Detector Reagent	101996	231312
SBG Reagent	102000	225108
Sample Diluent	102002	225109
Calibrators	103221	216401
Control 1	103222	216402
Control 2	103223	216402
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

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	216402	1.66	1.24-1.87
Control 2	216402	73.3	51.5-81.7
Panel 1	124607	3.09	2.54-3.81
Panel 2	208403	121	101-179
Panel 3	208403	2.22	1.81-3.42

Abeta 42

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	216402	4.20	3.18-4.77
Control 2	216402	94.8	83.0-127
Panel 1	124607	5.84	4.55-7.43
Panel 2	208403	7.36	6.37-10.7
Panel 3	208403	103	90.8-148

Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	216402	11.5	8.89-13.3
Control 2	216402	178	152-229
Panel 1	124607	132	114-171
Panel 2	208403	132	111-208
Panel 3	208403	83.5	67.4-122

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

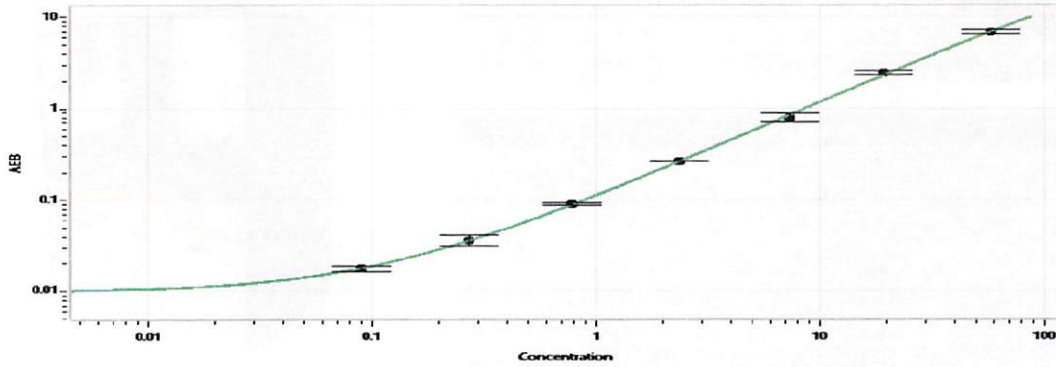
Review/Approval

Sonia Khillan
Name

QC
Title

Sonia Khillan 06DEC2022
Signature/Date

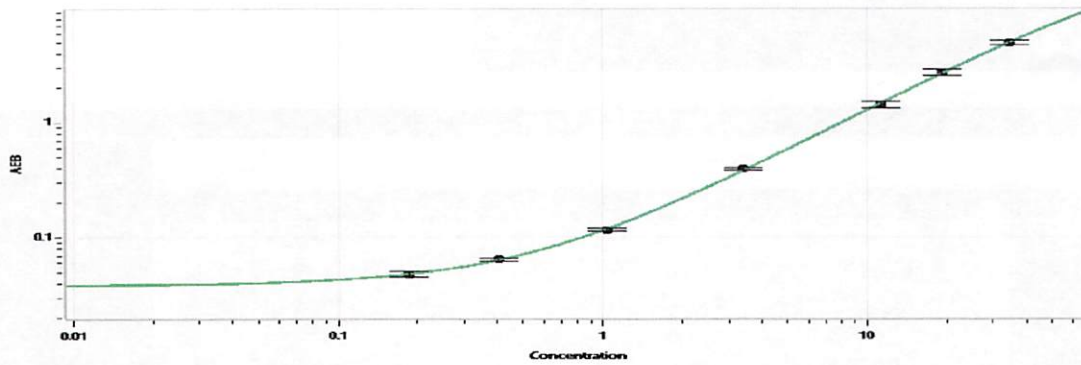
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.090
C	0.272
D	0.787
E	2.36
F	7.30
G	19.4
H	58.2

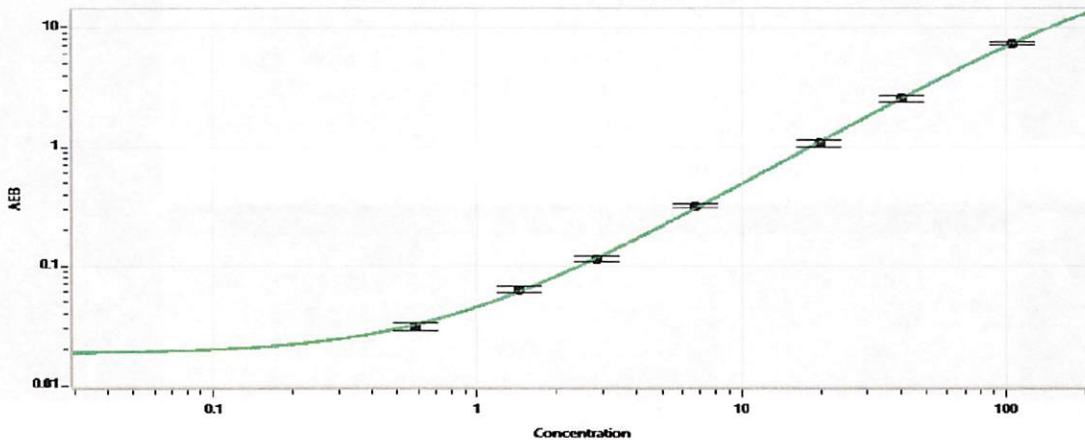
Abeta 42 Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.186
C	0.405
D	1.04
E	3.39
F	11.3
G	19.2
H	34.5

Abeta 40 Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.583
C	1.44
D	2.82
E	6.72
F	19.7
G	40.3
H	105