

Product Number:	101995
Lot Number:	503400
Expiration:	13-Apr-2023
Platform(s):	HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101998	210311
Detector Reagent	101996	214403
SBG Reagent	102000	212905
Sample Diluent	102002	212913
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	1.78	1.21-1.82
Control 2	216402	59.3	51.5-77.2
Panel 1	124607	3.36	2.54-3.81
Panel 2	124607	176	176-264
Panel 3	124607	5.15	3.61-5.86

Abeta 42

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	216402	3.91	3.12-4.68
Control 2	216402	93.7	86.7-130
Panel 1	124607	5.51	4.55-7.43
Panel 2	124607	5.69	4.54-7.13
Panel 3	124607	87.8	81.4-126

Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	216402	10.8	8.78-13.2
Control 2	216402	181	156-234
Panel 1	124607	131	114-171
Panel 2	124607	98.3	83.2-125
Panel 3	124607	169	136-209

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

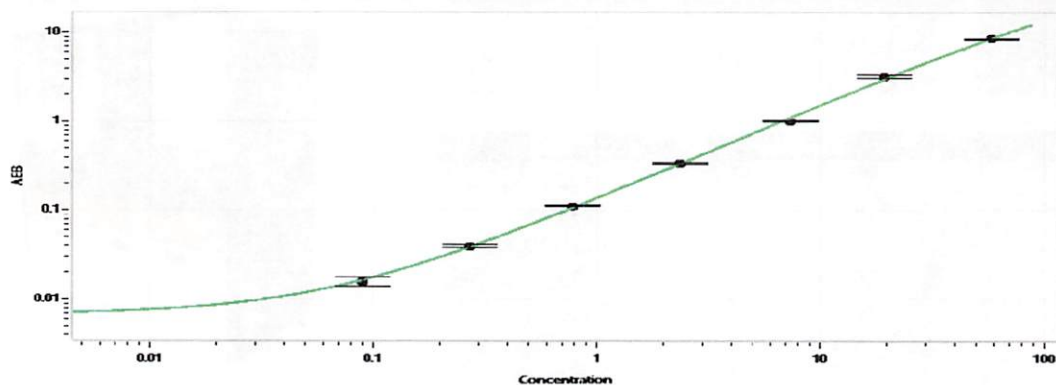
Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 08SEP2022
Signature/Date

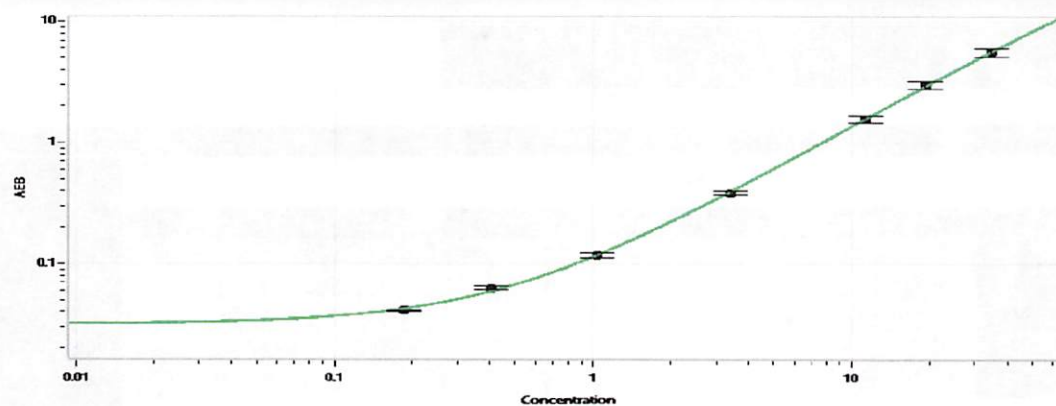
Tau Calibration Curve



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

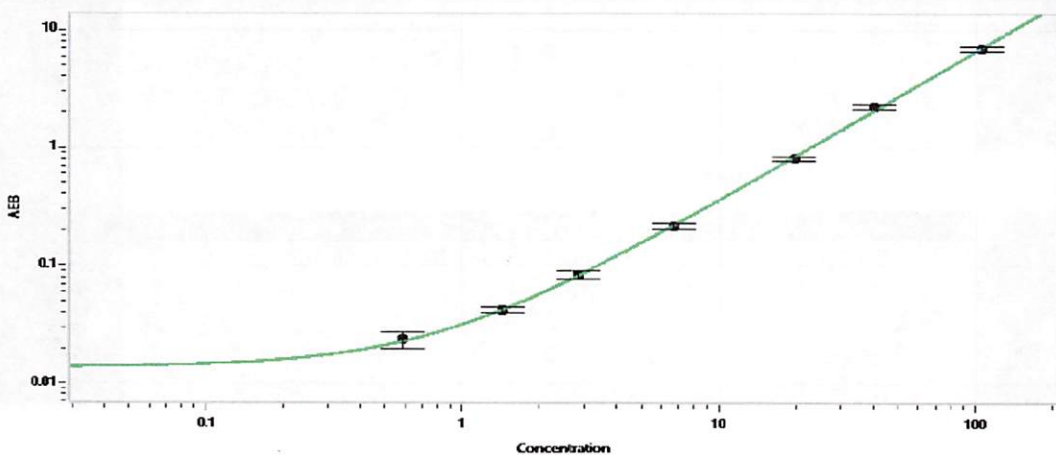
Abeta 42 Calibration Curve



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