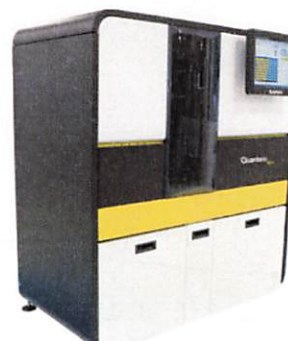


Product Number:	101995
Lot Number:	503705
Expiration:	7-Sep-2023
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101998	225009
Detector Reagent	101996	234806
SBG Reagent	102000	234906
Sample Diluent	102002	234907
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	1.88	1.17 - 2.07
Control 2	216402	61.4	50.9 - 80.2
Panel 1	211905	2.28	1.73 - 2.60
Panel 2	208403	122	96.9 - 175
Panel 3	208403	2.90	1.89 - 3.46

Abeta 42

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	216402	4.01	3.19 - 4.78
Control 2	216402	85.2	72.8 - 129
Panel 1	211905	1.91	1.32 - 2.42
Panel 2	208403	7.96	6.39 - 10.4
Panel 3	208403	99.8	82.7 - 148

Abeta 40

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	216402	11.9	9.01 - 13.5
Control 2	216402	151	135 - 230
Panel 1	211905	25.6	18.9 - 28.4
Panel 2	208403	126	98.1 - 208
Panel 3	208403	78.0	61.5 - 121

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

Review/Approval

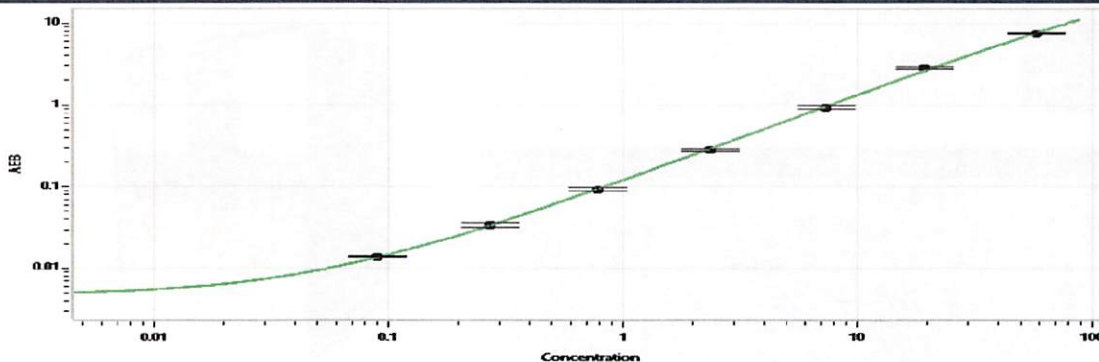
Sonia Khilkin
Name

QC

Title

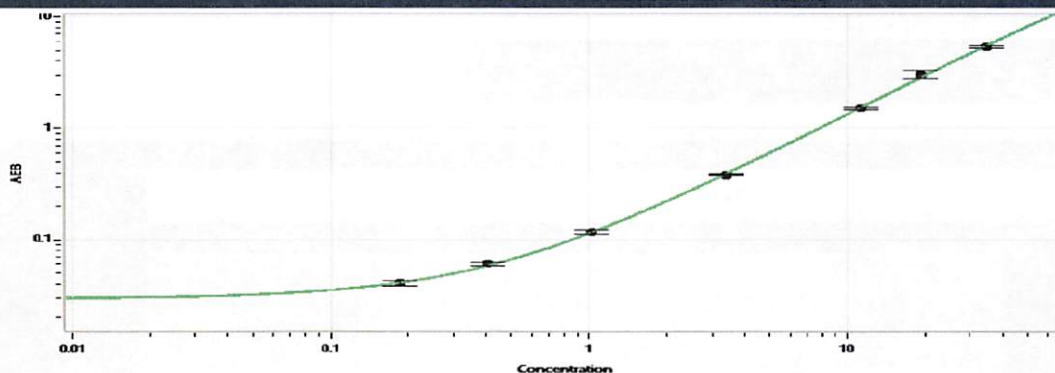
Sonia Khilkin / 21 DEC 2022
Signature/Date

Tau Calibration Curve



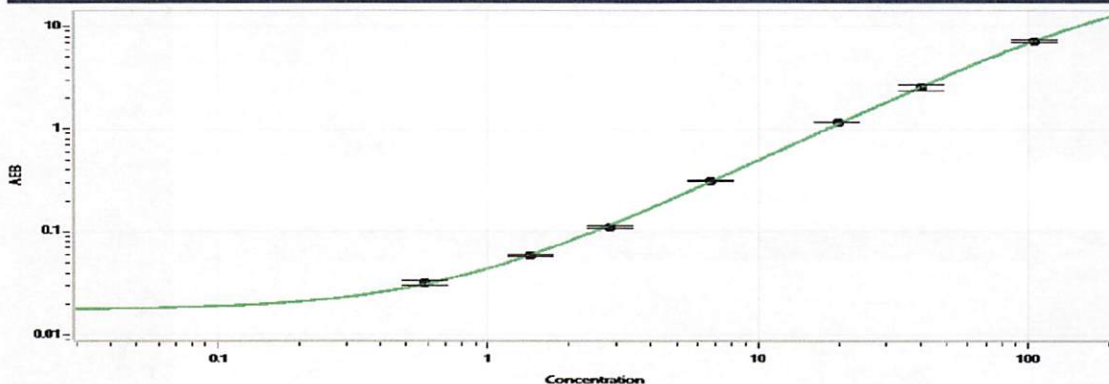
Calibrator Levels (pg/mL)	
A	0.000
B	0.090
C	0.272
D	0.787
E	2.36
F	7.30
G	19.4
H	58.2

Abeta 42 Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.186
C	0.405
D	1.03
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