

Product Number:	101995
Lot Number:	503793
Expiration:	2-Mar-2024
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	101998	306107	N/A
Detector Reagent	101996	306201	N/A
SBG Reagent	102000	306202	N/A
Sample Diluent	102002	306203	N/A
Calibrators	103221	317710	317710A
Control 1	103222	316407	316407A
Control 2	103223	316407	316407A
SBG Diluent (SRX)†	103158	N/A <sup>1</sup>	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>	N/A <sup>1</sup>

<sup>1</sup> 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)	pec. Range (pg/mL)
Control 1	316407A	1.53	1.24 - 1.85
Control 2	316407A	63.2	53.2 - 79.7
Panel 1	124607	2.67	2.54 - 3.81
Panel 2	208403	126	86.2 - 177
Panel 3	208403	2.15	1.91 - 3.59

**Abeta 42**

	Lot	Result (pg/mL)	pec. Range (pg/mL)
Control 1	316407A	2.07	1.59 - 2.44
Control 2	316407A	80.5	60.8 - 91.1
Panel 1	124607	5.22	4.55 - 7.43
Panel 2	208403	7.18	6.55 - 10.3
Panel 3	208403	94.0	80.8 - 145

**Abeta 40**

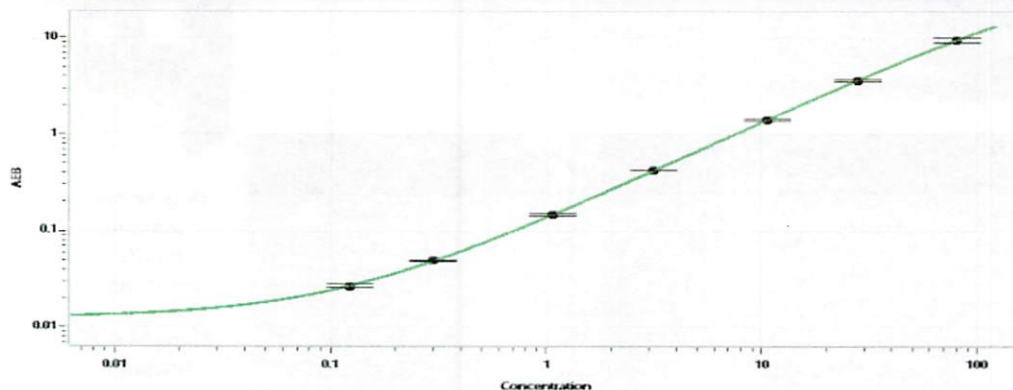
	Lot	Result (pg/mL)	pec. Range (pg/mL)
Control 1	316407A	5.42	3.87 - 6.08
Control 2	316407A	130	109 - 164
Panel 1	124607	123	114 - 171
Panel 2	208403	128	101 - 203
Panel 3	208403	78.7	63.8 - 118

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

Review/Approval

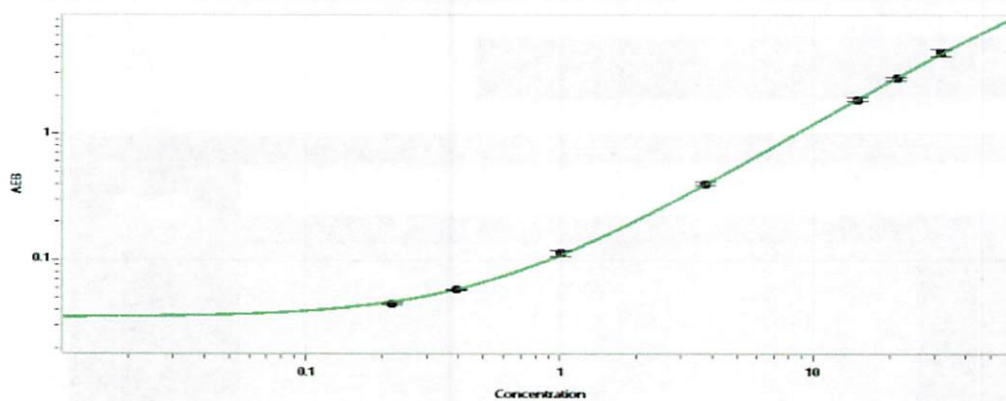
Sonia Khilkin QC Sonia 12/1/23 18 JUL 2023  
Name Title Signature/Date

**Tau Calibration Curve**



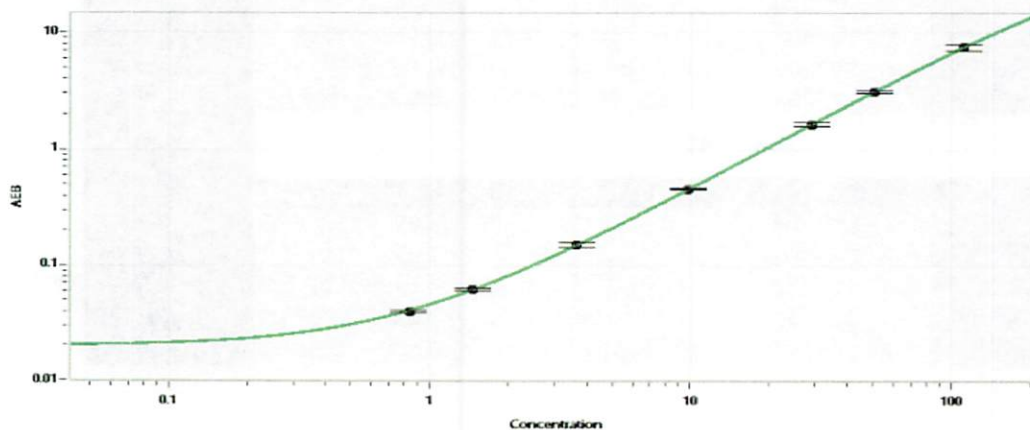
Calibrator Levels (pg/mL)	
A	0.000
B	0.123
C	0.300
D	1.07
E	3.16
F	10.5
G	27.8
H	80.0

**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