

Product Number:	101995
Lot Number:	503239
Expiration:	21-Aug-2022
Platform(s):	HD-1/HD-X & SR-X



Data below represents results generated on the Simoa® HD-1 / HD-X / SR-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101998	133101
Detector Reagent	101996	134209
SBG Reagent	102000	133206
Sample Diluent	102002	123311
Calibrators	103221	134903
Control 1	103222	134904
Control 2	103223	134904
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)	Mean Range (pg/mL)
Control 1	134904	2.91	2.33-3.49
Control 2	134904	112	98.2-147
Panel 1	211905	2.25	1.70-2.55
Panel 2	124607	217	176-264
Panel 3	124607	5.42	3.61-5.86

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134904	5.25	4.13-6.98
Control 2	134904	138	121-186
Panel 1	211905	2.00	1.27-2.34
Panel 2	124607	5.65	4.54-7.13
Panel 3	124607	103	81.4-126

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134904	14.2	12.3-18.4
Control 2	134904	231	204-330
Panel 1	211905	22.0	18.2-27.3
Panel 2	124607	106	83.2-125
Panel 3	124607	167	136-209

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

Review/Approval

Matt Smart

QC Director

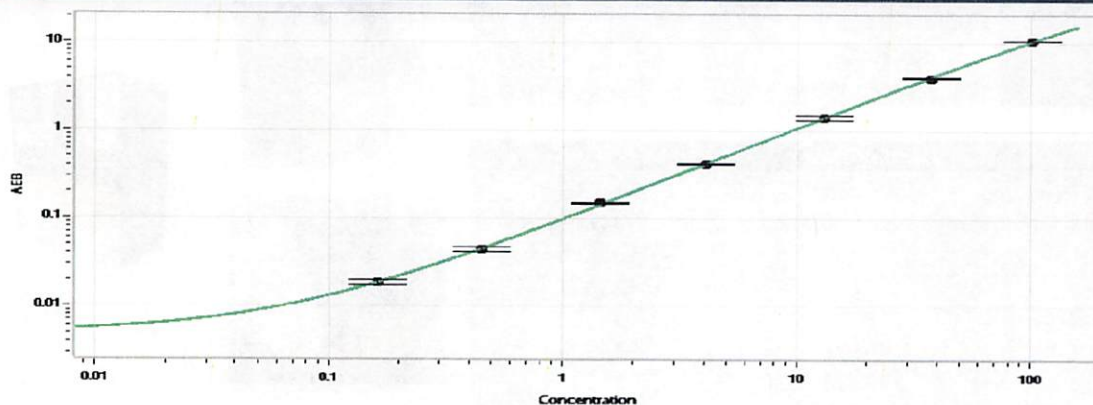
[Handwritten Signature] 06 Jun 2022

Name

Title

Signature/Date

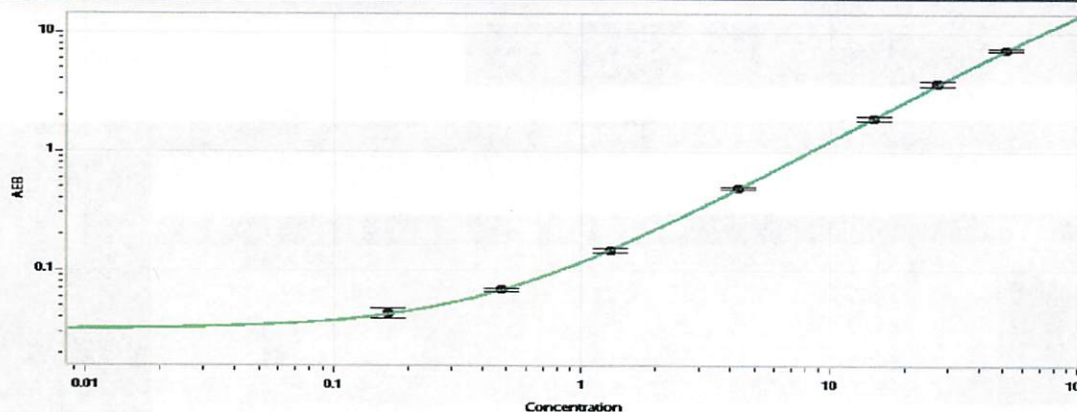
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.162
C	0.455
D	1.44
E	4.07
F	13.0
G	37.5
H	101

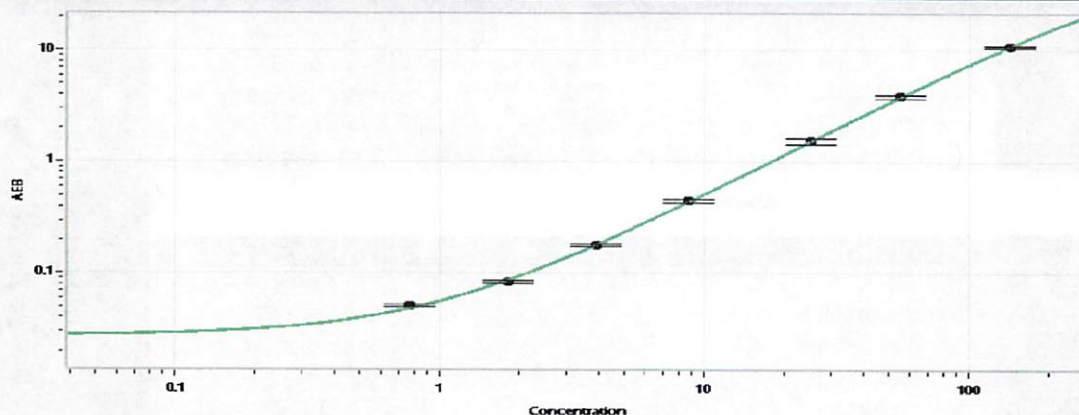
Abeta 42



Calibrator Levels (pg/mL)

A	0.000
B	0.166
C	0.474
D	1.30
E	4.25
F	15.1
G	27.3
H	51.6

Abeta 40



Calibrator Levels (pg/mL)

A	0.000
B	0.774
C	1.81
D	3.89
E	8.71
F	25.4
G	55.2
H	142