

Product Number:	101995
Lot Number:	503519
Expiration:	10-Nov-2022
Platform(s):	HD-1/HD-X & SR-X



Component	Part Number	Lot Number
Bead Reagent	101998	133101
Detector Reagent	101996	134209
SBG Reagent	102000	133206
Sample Diluent	102002	216202
Calibrators	103221	134903
Control 1	103222	134904
Control 2	103223	134904
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

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

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	2.63	2.28-3.42
Control 2	134904	109	94.0-146
Panel 1	211905	2.21	1.72-2.57
Panel 2	124607	194	176-264
Panel 3	124607	5.33	3.61-5.86

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134904	5.17	4.12-6.84
Control 2	134904	136	115-184
Panel 1	211905	2.08	1.30-2.43
Panel 2	124607	5.84	4.54-7.13
Panel 3	124607	107	81.4-126

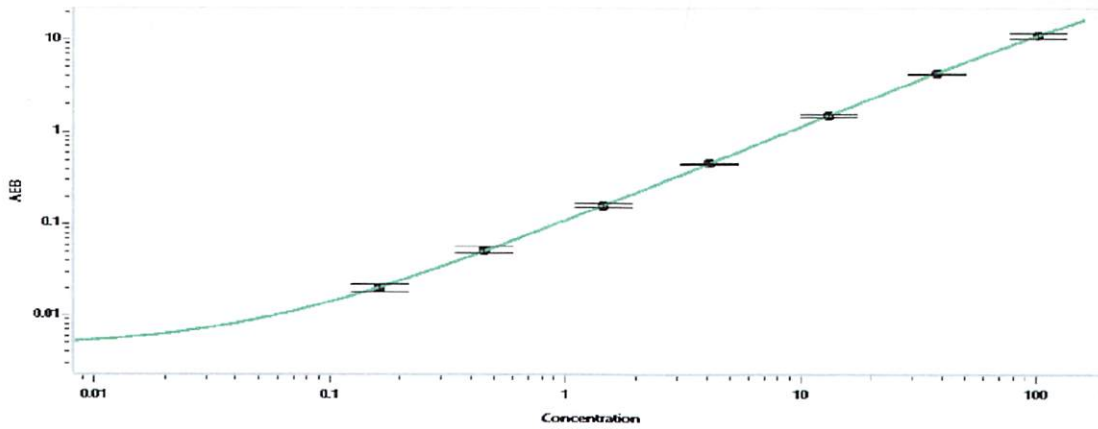
Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134904	14.5	12.2-18.2
Control 2	134904	251	205-323
Panel 1	211905	25	18.6-28.0
Panel 2	124607	103	83.2-125
Panel 3	124607	178	136-209

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

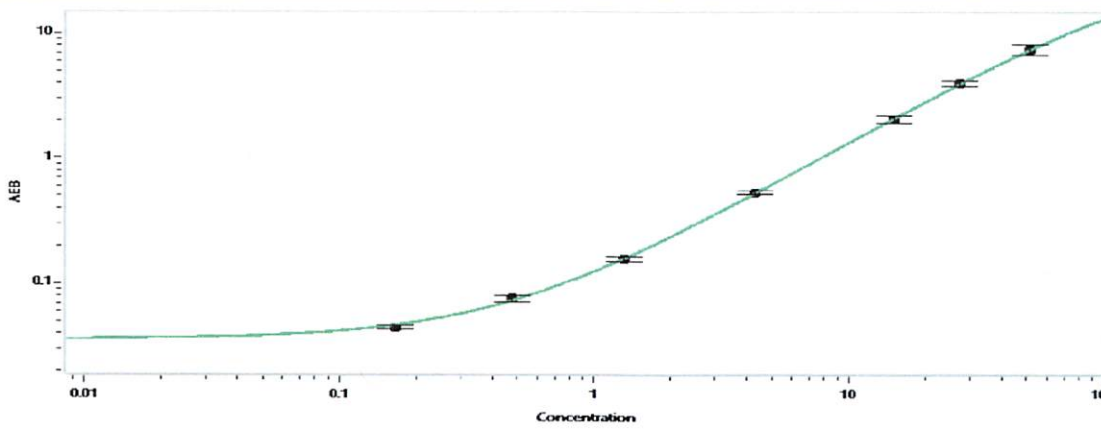
Review/Approval	Matt Smart	QC Director		24 Jun 2022
	Name	Title	Signature/Date	

Tau Calibration Curve



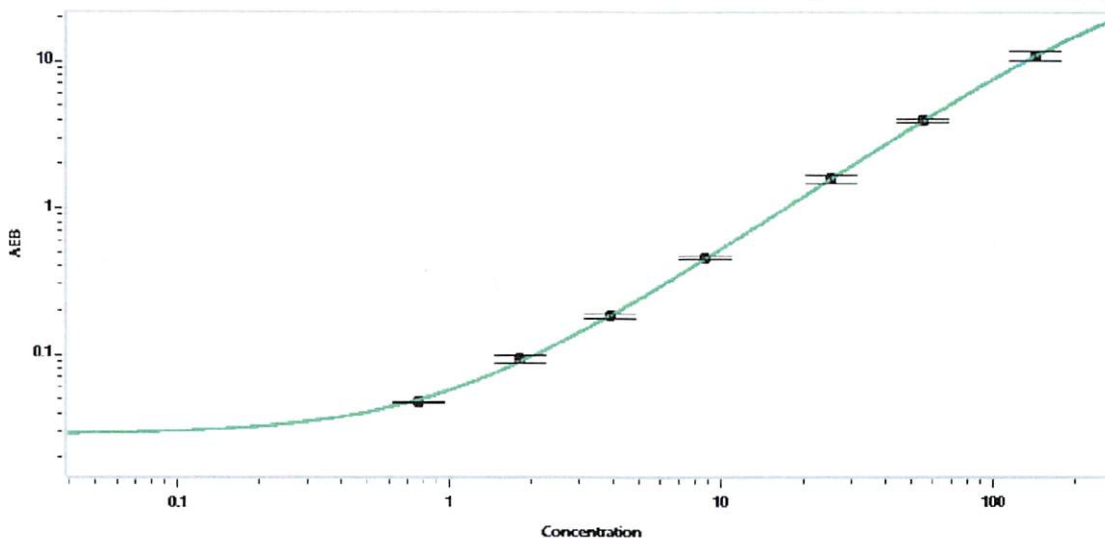
Calibrator Levels (pg/mL)	
A	0.000
B	0.162
C	0.455
D	1.44
E	4.07
F	13.0
G	37.5
H	101

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