

Product Number:	101995
Lot Number:	503245
Expiration:	5-Aug-2022
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	129125
Detector Reagent	101996	128904
SBG Reagent	102000	128905
Sample Diluent	102002	123412
Calibrators	103221	128302
Control 1	103222	128303
Control 2	103223	128303
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	128303	2.97	2.55 - 4.07
Control 2	128303	113	97.6 - 146
Panel 1	124607	3.16	2.57 - 3.86
Panel 2	124607	221	175 - 263
Panel 3	124607	4.70	3.83 - 5.76

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	128303	5.01	4.27 - 6.40
Control 2	128303	123	110 - 166
Panel 1	124607	5.58	4.44 - 7.68
Panel 2	124607	5.63	4.75 - 7.24
Panel 3	124607	92.0	85.2 - 128

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	128303	14.3	11.6 - 17.5
Control 2	128303	209	149 - 351
Panel 1	124607	131	116 - 174
Panel 2	124607	96.6	84.9 - 127
Panel 3	124607	150	140 - 213

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

Review/Approval

Shivani Goel

QC Manager

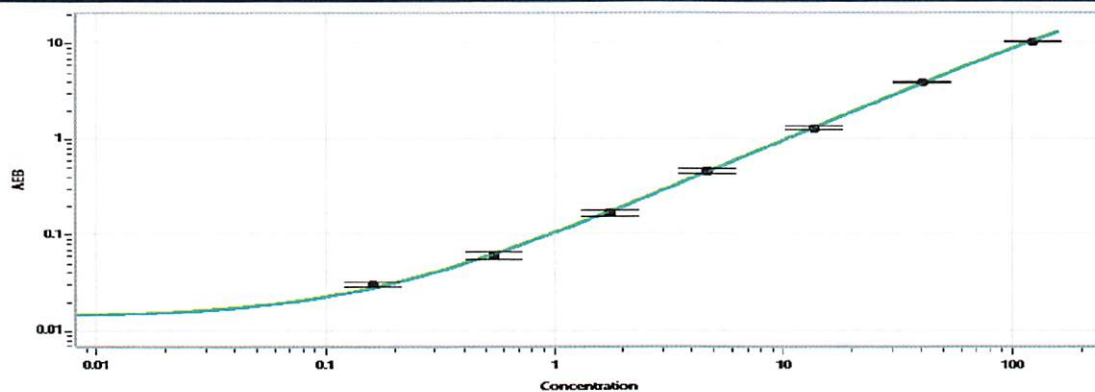
Shwam Goel 02 May 2022

Name

Title

Signature/Date

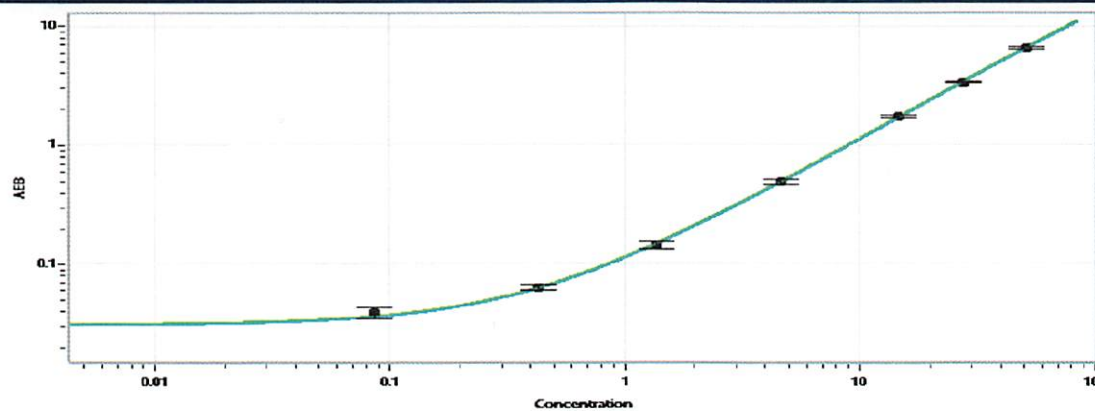
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.161
C	0.546
D	1.75
E	4.65
F	13.6
G	40.4
H	123

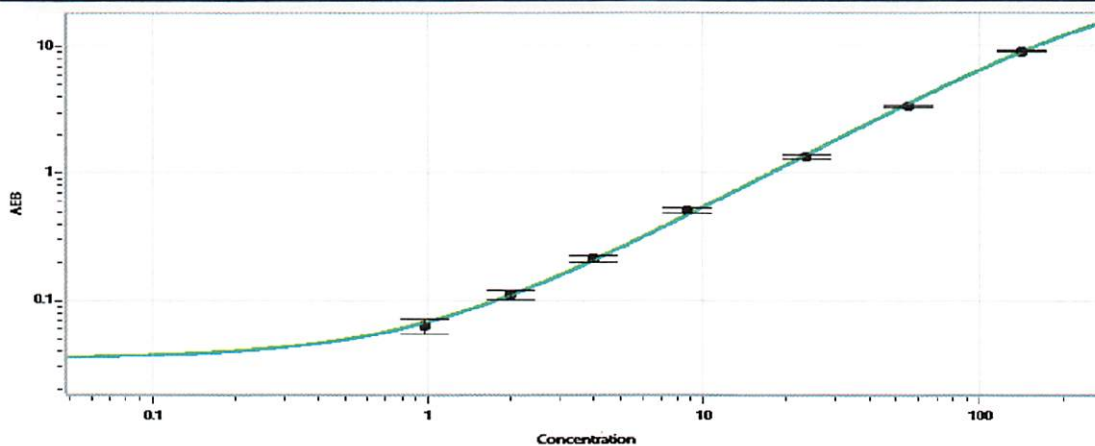
Abeta 42



Calibrator Levels (pg/mL)

A	0.000
B	0.086
C	0.428
D	1.37
E	4.65
F	14.8
G	27.7
H	51.2

Abeta 40



Calibrator Levels (pg/mL)

A	0.000
B	0.966
C	1.98
D	3.99
E	8.72
F	23.5
G	55.0
H	143