

Product Number:	101995
Lot Number:	502238
Expiration:	21-Nov-2020
Platform(s):	HD-1/HD-X & SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101998	001513
Detector Reagent	101996	935301
SBG Reagent	102000	934017
Sample Diluent	102002	934016
Calibrators	103221	003701 & 004301
Controls	103222/103223	003702
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	003702	2.11	1.39-4.16
Control 2	003702	89.9	50.0-150
Panel 1	003802	1.39	0.889-2.57
Panel 2	003802	133	105-194
Panel 3	003802	0.730	0.498-1.55

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	003702	4.34	2.77-6.94
Control 2	003702	117	81.3-210
Panel 1	003802	4.33	1.27-8.14
Panel 2	003802	3.92	1.24-7.93
Panel 3	003802	97.9	73.9-169

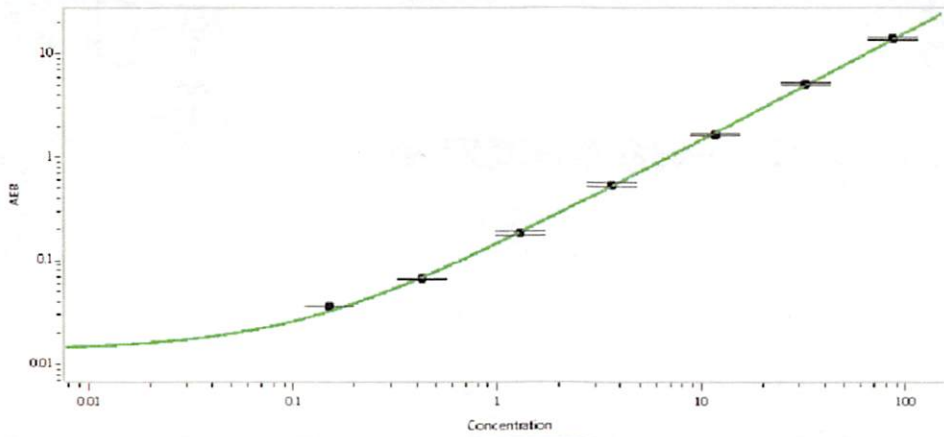
Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	003702	15.0	9.40-21.9
Control 2	003702	231	149-347
Panel 1	003802	62.1	44.3-107
Panel 2	003802	31.9	25.5-52.5
Panel 3	003802	70.3	45.2-132

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

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>p. stadler</i> 24 Feb 2020
	Name	Title	Signature/Date

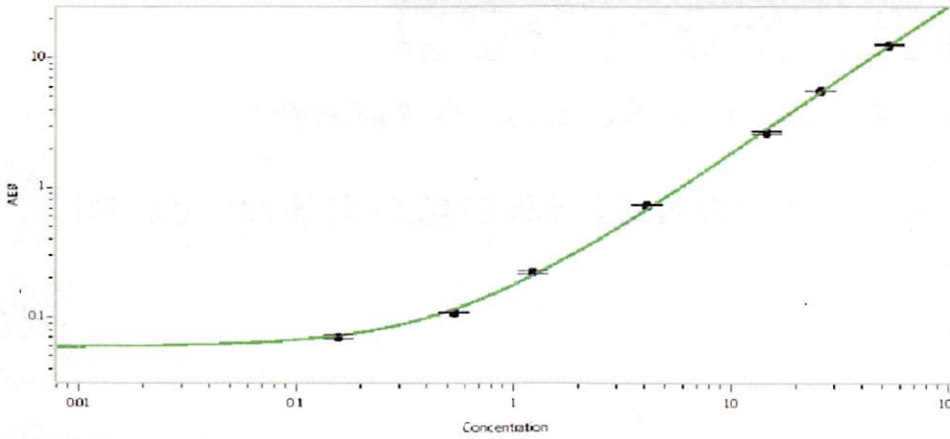
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.151
C	0.426
D	1.30
E	3.65
F	11.7
G	32.2
H	87.0

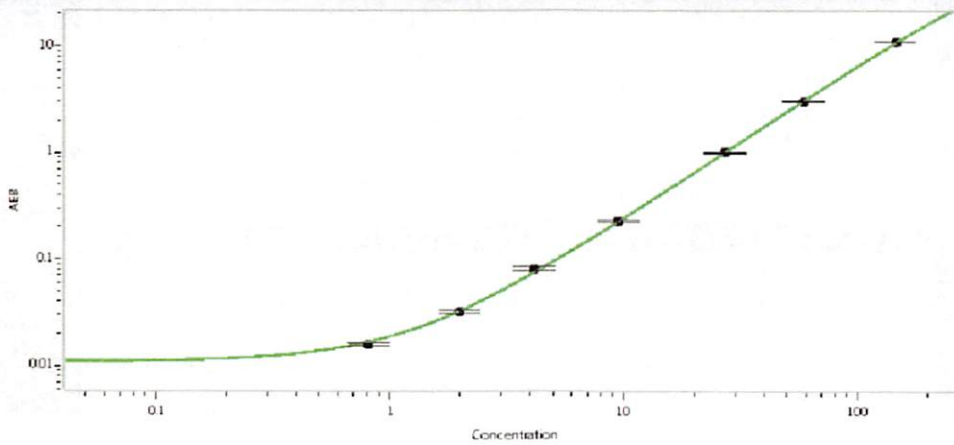
Abeta 42



Calibrator Levels (pg/mL)

A	0.000
B	0.158
C	0.538
D	1.24
E	4.14
F	14.6
G	25.9
H	53.7

Abeta 40



Calibrator Levels (pg/mL)

A	0.000
B	0.808
C	1.98
D	4.16
E	9.46
F	26.9
G	58.9
H	145