

Product Number:	101995
Lot Number:	502593
Expiration:	23-Jun-2021
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101998	029713
Detector Reagent	101996	029714
SBG Reagent	102000	030005
Sample Diluent	102002	030003
Calibrators	103221	027402
Control 1	103222	931002
Control 2	103223	931002
SBG Diluent (SRX)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	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).



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

**Release Materials**

**Tau**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931002	1.80	0.775-2.26
Control 2	931002	82.0	40.8-92.3
Panel 1	003802	1.52	0.935-2.42
Panel 2	003802	161	106-194
Panel 3	003802	1.29	0.552-1.68

**Abeta 42**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931002	5.17	4.47-6.79
Control 2	931002	132	107-160
Panel 1	003802	4.51	2.66-6.02
Panel 2	003802	4.27	2.78-5.20
Panel 3	003802	121	88.7-140

**Abeta 40**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931002	19.5	13.6-20.9
Control 2	931002	243	203-305
Panel 1	003802	75.8	47.2-101
Panel 2	003802	41.3	26.4-50.9
Panel 3	003802	101	50.4-128

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

Review/Approval

Shivani Goel

QC Manager

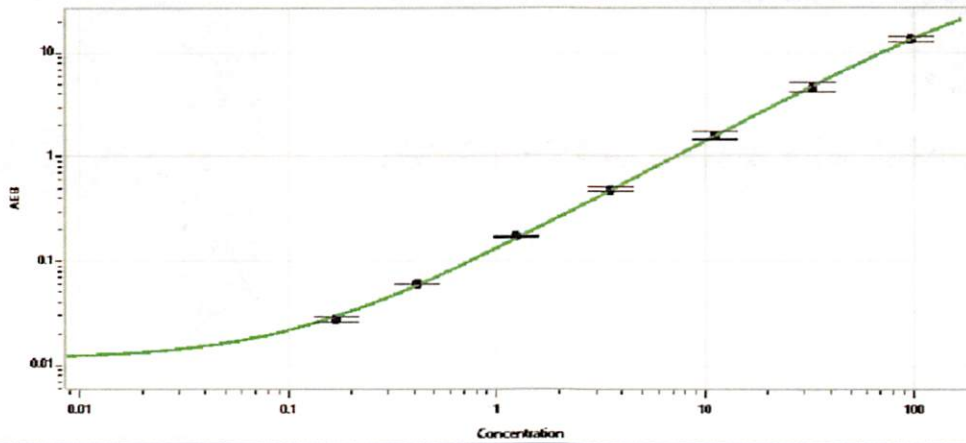
Shivani Goel 11 Nov 2020

Name

Title

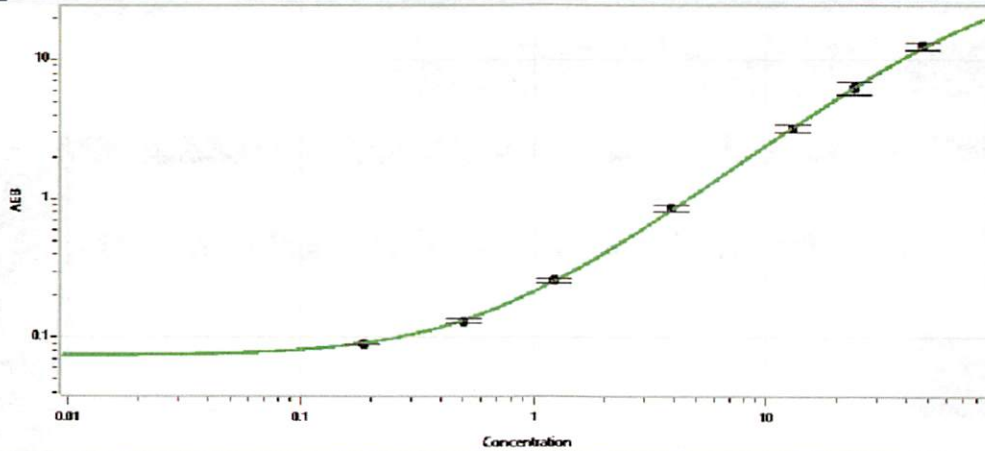
Signature/Date

**Tau Calibration Curve**



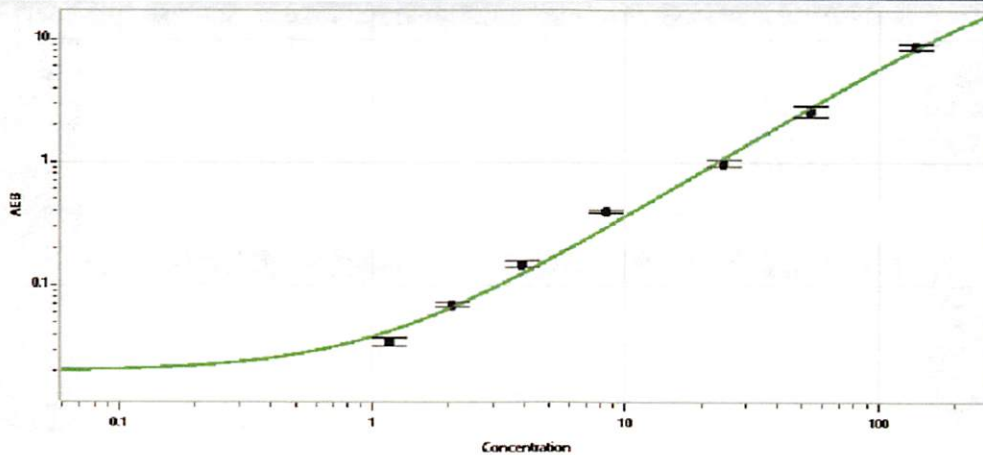
Calibrator Levels (pg/mL)	
A	0.000
B	0.170
C	0.413
D	1.24
E	3.52
F	11.2
G	32.7
H	98.2

**Abeta 42**



Calibrator Levels (pg/mL)	
A	0.000
B	0.186
C	0.501
D	1.23
E	3.91
F	13.0
G	24.0
H	47.3

**Abeta 40**



Calibrator Levels (pg/mL)	
A	0.000
B	1.17
C	2.08
D	3.93
E	8.36
F	24.5
G	54.0
H	141