

Product Number:	101995
Lot Number:	502701
Expiration:	11-Jan-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	101909
Detector Reagent	101996	105002
SBG Reagent	102000	102009
Sample Diluent	102002	102011
Calibrators	103221	102503
Control 1	103222	102504
Control 2	103223	102504
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).

**Release Materials**

**Tau**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	102504	1.74	1.29-1.94
Control 2	102504	67.1	46.2-85.9
Panel 1	003802	2.30	0.898-2.55
Panel 2	003802	154	104-190
Panel 3	003802	1.56	0.809-1.54

**Abeta 42**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	102504	5.04	3.56-5.66
Control 2	102504	118	91.9-148
Panel 1	003802	4.77	2.54-6.08
Panel 2	003802	4.53	2.67-5.20
Panel 3	003802	103	88.4-138

**Abeta 40**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	102504	14.3	8.74-17.8
Control 2	102504	201	160-246
Panel 1	003802	68.9	44.5-103
Panel 2	003802	36.4	25.1-51.1
Panel 3	003802	80.2	46.9-127

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

Review/Approval

Shivani Goel

QC Manager

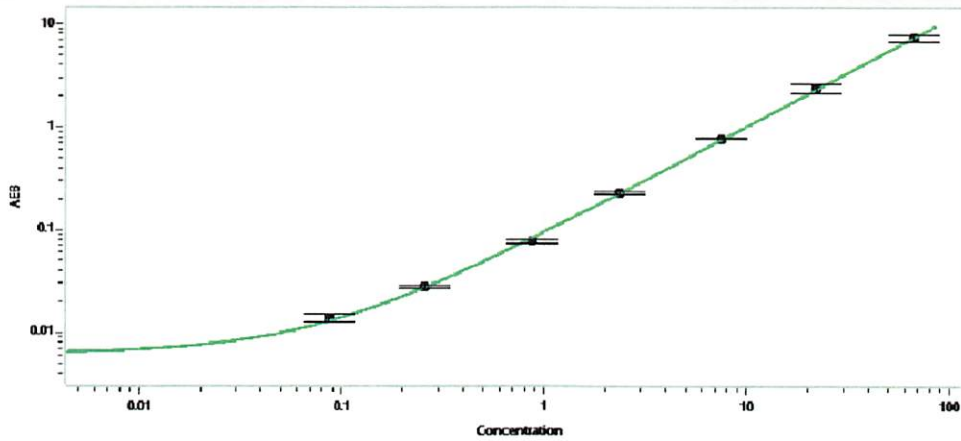
*Shivani Goel 12 May 2022*

Name

Title

Signature/Date

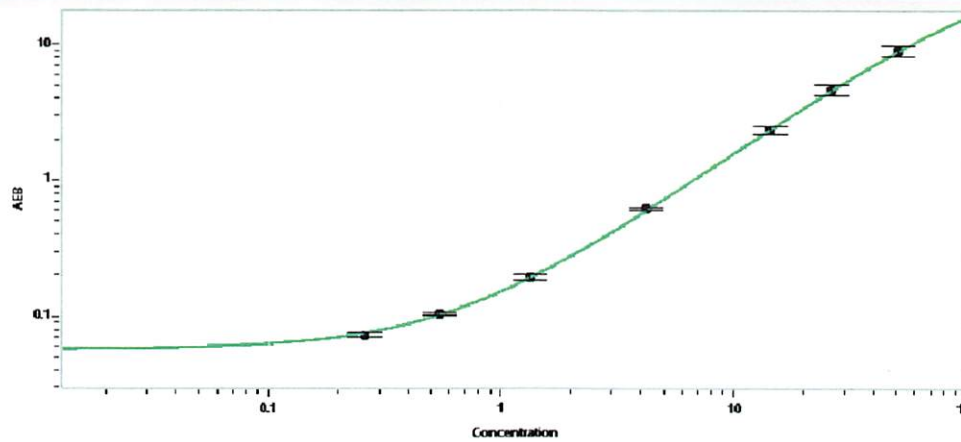
**Tau Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.086
C	0.255
D	0.867
E	2.35
F	7.52
G	22.0
H	66.0

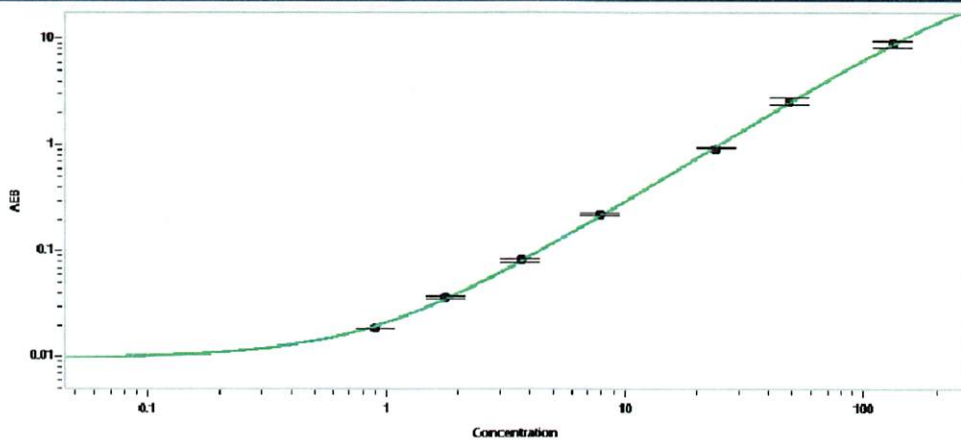
**Abeta 42**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.257
C	0.545
D	1.32
E	4.22
F	14.3
G	26.3
H	50.6

**Abeta 40**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.900
C	1.78
D	3.66
E	7.88
F	23.9
G	48.7
H	132