

Product Number:	101995
Lot Number:	501643
Expiration:	22-Jan-2020
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101998	903807
Detector Reagent	101996	903901
SBG Reagent	102000	902802
Sample Diluent	102002	902803
Calibrators	103221	820001
Controls	103222-103223	834609
SBG Diluent (SR-X)†	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 SR-X Kit Instructions (KI-0083).

Release Materials

Tau

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	834609	2.21	1.60-2.41
Control 2	834609	88.4	67.2-101
Panel 1	MF-062018-04	1.31	1.16-2.04
Panel 2	MF-010818-02	181	156-234
Panel 3	MF-010818-03	3.64	1.30-5.98

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	834609	5.00	3.33-6.56
Control 2	834609	113	87.7-132
Panel 1	MF-062018-04	6.60	5.28-7.92
Panel 2	MF-010818-02	8.89	7.07-10.7
Panel 3	MF-010818-03	119	96.1-150

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	834609	17.0	10.2-18.1
Control 2	834609	230	172-259
Panel 1	MF-062018-04	116	102-158
Panel 2	MF-010818-02	133	107-200
Panel 3	MF-010818-03	87.2	71.5-124

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

Review/Approval

Patrizia Stadler Assoc. Director, QC

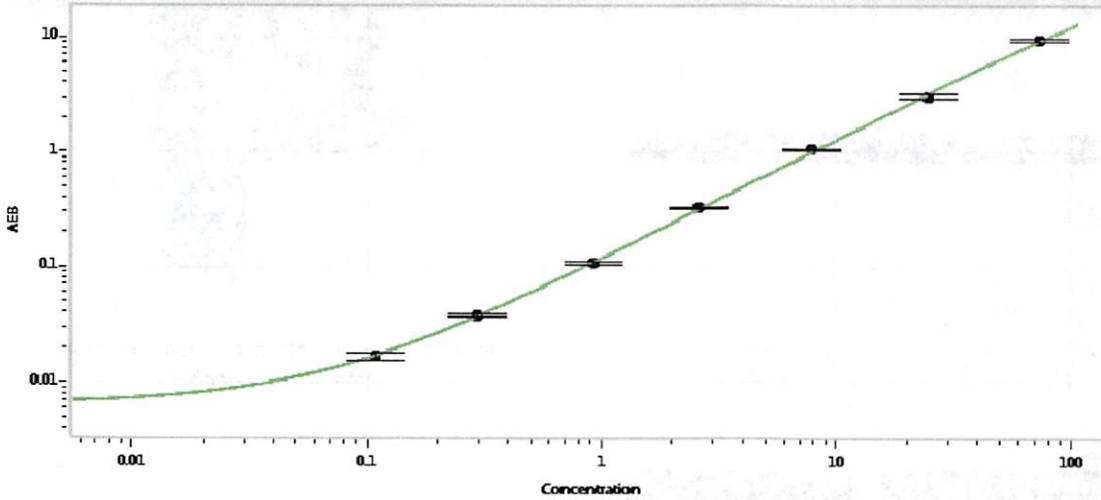
P. Sooler 11 Feb 2019

Name Title Signature/Date

Simoa™ Neuro 3-Plex A Kit

Certificate of Analysis

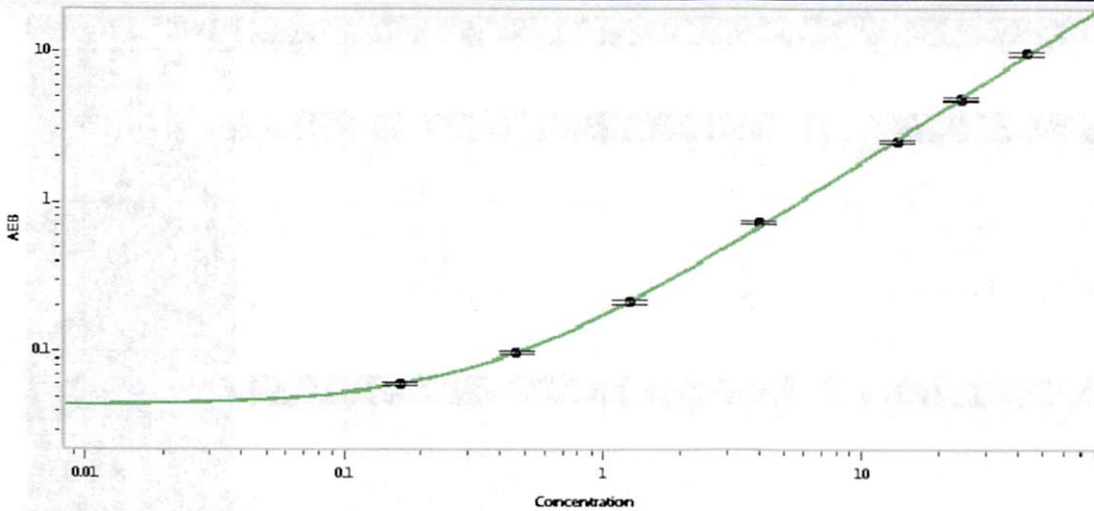
Tau Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.109
C	0.294
D	0.919
E	2.59
F	7.88
G	24.7
H	73.5

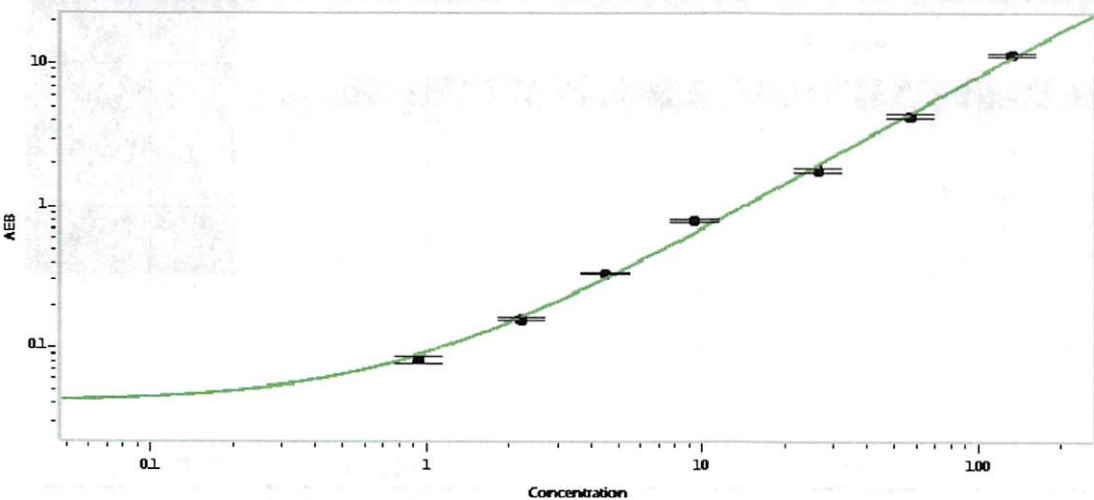
Abeta 42 Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.165
C	0.462
D	1.27
E	4.03
F	13.7
G	24.2
H	43.1

Abeta 40 Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.938
C	2.22
D	4.48
E	9.43
F	26.3
G	57.3
H	133