

Product Number:	101995
Lot Number:	501965
Expiration:	13-Jun-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	921704
Detector Reagent	101996	921708
SBG Reagent (HD-1)	102000	921203
Sample Diluent	102002	921107
Calibrators	103221	915601 & 916501
Controls	103222/103223	915602
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	915602	1.65	1.32-1.99
Control 2	915602	71.7	55.7-83.5
Panel 1	MF-062018-04	1.29	1.18-2.03
Panel 2	MF-010818-02	171	155-233
Panel 3	MF-010818-03	3.69	1.30-5.98

Abeta 42

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	915602	5.26	4.09-6.13
Control 2	915602	123	94.4-148
Panel 1	MF-062018-04	7.17	5.32-7.98
Panel 2	MF-010818-02	10.6	6.80-11.4
Panel 3	MF-010818-03	135	96.1-150

Abeta 40

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	915602	13.6	11.0-16.5
Control 2	915602	231	184-277
Panel 1	MF-062018-04	143	104-159
Panel 2	MF-010818-02	148	122-192
Panel 3	MF-010818-03	105	71.5-124

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

Review/Approval	Daniel Ritchie	Senior QC Analyst	<i>[Signature]</i> 25 Aug 2018
	Name	Title	Signature/Date

*Signing on behalf of Patrizia Stadler