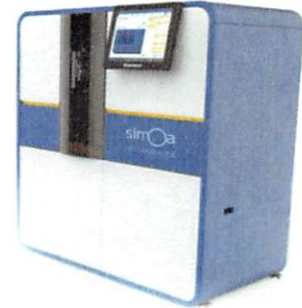


|                 |                  |
|-----------------|------------------|
| Product Number: | 101552           |
| Lot Number:     | 501939           |
| Expiration:     | 23-Mar-2020      |
| Platform(s):    | HD-1/HD-x & SR-X |



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

| Component            | Part Number   | Lot Number       |
|----------------------|---------------|------------------|
| Bead Reagent         | 101554        | 908805           |
| Detector Reagent     | 101553        | 908509           |
| SBG Reagent          | 101555        | 908508           |
| Sample Diluent       | 101556        | 908204           |
| Calibrators          | 101559        | 912902           |
| Controls             | 101560-101561 | 912903           |
| SBG Diluent (SR-X) † | 103158        | N/A <sup>1</sup> |
| RGP Reagent          | 103159        | N/A <sup>1</sup> |

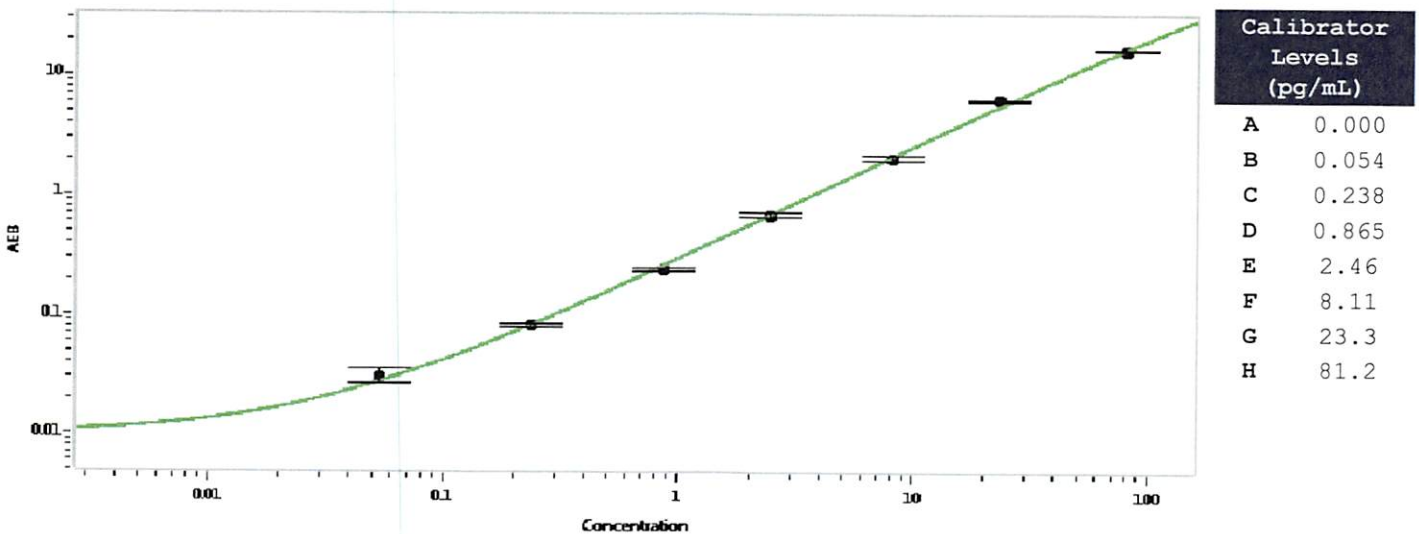
<sup>1</sup> RGP is not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ Human Total Tau Advantage SR-X Kit Instructions (KI-0034).

| Release Materials* |              |                |                    |
|--------------------|--------------|----------------|--------------------|
|                    | Lot          | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1          | 912903       | 0.766          | 0.680-1.28         |
| Control 2          | 912903       | 53.5           | 42.5-63.8          |
| Panel 1            | MF-013017-02 | 0.942          | 0.251-1.23         |
| Panel 2            | MF-050317-13 | 1.36           | 0.326-2.04         |
| Panel 3            | MF-050417-01 | 69.3           | 38.2-115           |
| Panel 4            | MF-050417-02 | 165            | 99.8-299           |

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

Calibration Curve



Review/Approval

Daniel Ritchie Senior QC Analyst

Name

Title

*[Signature]* / 22 Aug 2019  
Signature/Date

\*Signing on behalf of Patrizia Stadler