

|                         |                    |
|-------------------------|--------------------|
| <b>Product Number:</b>  | 101552             |
| <b>Lot Number:</b>      | 504003             |
| <b>Expiration Date:</b> | 08-Nov-2024        |
| <b>Platform(s):</b>     | SR-X and HD-1/HD-X |



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

| Component           | Part Number | Lot Number       |
|---------------------|-------------|------------------|
| Bead Reagent        | 101554      | 404711           |
| Detector Reagent    | 101553      | 401816           |
| SBG Reagent         | 101555      | 401817           |
| Sample Diluent      | 101556      | 401905           |
| Calibrators         | 101559      | 403004           |
| Control 1           | 101560      | 403005           |
| Control 2           | 101561      | 403005           |
| SBG Diluent (SR-X)† | 103158      | N/A <sup>1</sup> |
| RGP Reagent         | 103159      | N/A <sup>1</sup> |

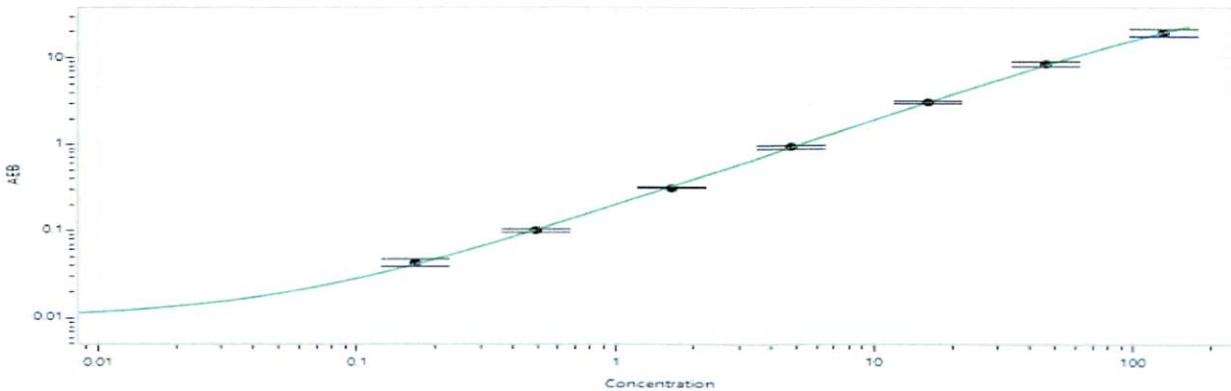
<sup>1</sup> Reagents are not Kit Lot Specific

† Specific to use on the SRX platform.  
Refer to the Tau Advantage SR-X Kit Instructions (KI-0034).

| Release Materials* |        |                |                     |
|--------------------|--------|----------------|---------------------|
|                    | Lot    | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1          | 403005 | 1.44           | 1.26 - 1.89         |
| Control 2          | 403005 | 83.5           | 71.7 - 108          |
| Panel 1            | 232607 | 0.408          | 0.214 - 1.16        |
| Panel 2            | 232607 | 0.229          | 0.014 - 1.16        |
| Panel 3            | 234908 | 18.2           | 14.7 - 22.1         |
| Panel 4            | 234908 | 30.9           | 22.3 - 33.4         |

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

**Calibration Curve**



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 0.167 |
| C                         | 0.493 |
| D                         | 1.64  |
| E                         | 4.78  |
| F                         | 16.1  |
| G                         | 46.0  |
| H                         | 130   |

Review/Approval

Cody Carlson

QC Manager

*[Signature]* 22 FEB 2024

Name

Title

Signature/Date