

<b>Product Number:</b>	101552
<b>Lot Number:</b>	504283
<b>Expiration Date:</b>	18Jul2025
<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

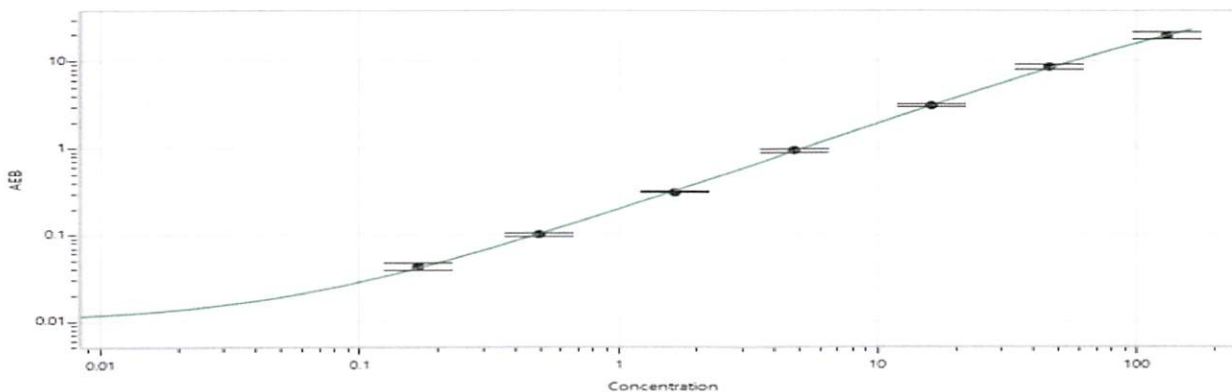
† Specific to use on the SRX platform.

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

Name

Title

Signature/Date

*[Handwritten Signature]* 05Jun2024