

<b>Product Number:</b>	101552
<b>Lot Number:</b>	504227
<b>Expiration Date:</b>	12-Aug-25
<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	411301
Detector Reagent	101553	407345
SBG Reagent	101555	411302
Sample Diluent	101556	404320
Calibrators	101559	416509
Control 1	101560	411725
Control 2	101561	411725
SBG Diluent (SRX)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

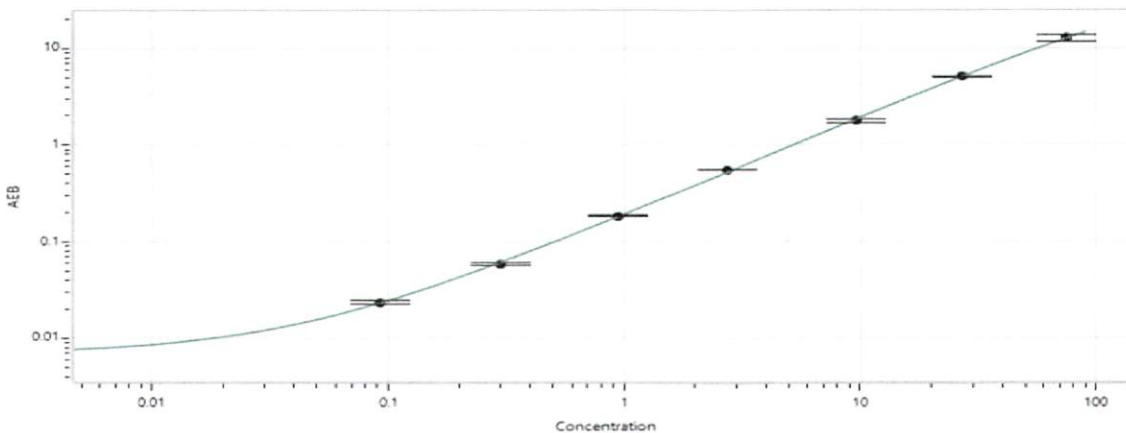
<sup>1</sup> Reagents are not Kit Lot Specific

† 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	411725	1.82	1.35 - 2.23
Control 2	411725	90.8	69.7 - 126
Panel 1	232607	0.480	0.172 - 1.12
Panel 2	232607	0.416	0.011 - 1.10
Panel 3	234908	18.5	14.7 - 22.1
Panel 4	234908	29.2	22.5 - 33.7

\* 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	0.092
C	0.301
D	0.937
E	2.73
F	9.60
G	27.1
H	74.7

Review/Approval

Name

Title

Signature/Date

25 JUN 2024