

Product Number:	101552
Lot Number:	503945
Expiration Date:	1-Aug-24
Platform(s):	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	332004
Detector Reagent	101553	321312
SBG Reagent	101555	325902
Sample Diluent	101556	325903
Calibrators	101559	325501
Control 1	101560	325502
Control 2	101561	325502
SBG Diluent (SRX)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

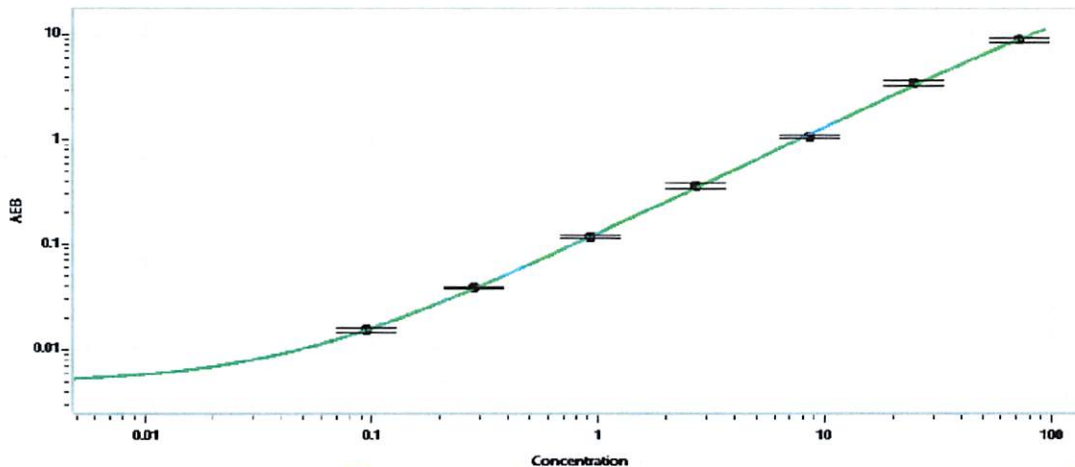
<sup>1</sup> Reagents are not Kit Lot Specific

† 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	325502	0.949	0.658 - 1.06
Control 2	325502	48.4	46.0 - 69.1
Panel 1	232607	0.695	0.514 - 1.05
Panel 2	232607	0.613	0.454 - 0.962
Panel 3	234908	16.9	14.8 - 22.1
Panel 4	234908	25.7	21.5 - 32.2

\* 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	0.095
C	0.283
D	0.928
E	2.68
F	8.58
G	24.7
H	71.6

Review/Approval

ROBERT SOMOGYI

QC MANAGER

RA [Signature] 18DEC2023

Name

Title

Signature/Date