

Product Number:	101552
Lot Number:	504322
Expiration Date:	1-Feb-25
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 Reagent	101556	325903
Reference Calibrators	101559	325501
Control 1	101560	325502
Control 2	101561	325502
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

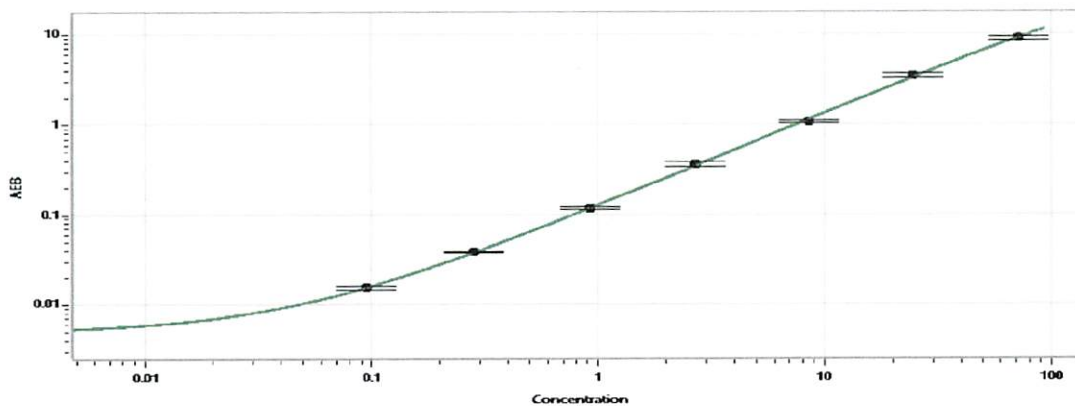
† Specific to use on the SR-X platform.

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

* 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

Cody Carlson

QC Manager

[Signature] 26 Jun 2024

Name

Title

Signature/Date