

Product Number:	101552
Lot Number:	504301
Expiration Date:	1-Oct-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	417904
Detector Reagent	101553	417802
SBG Reagent	101555	417905
Sample Diluent Reagent	101556	409215
Reference Calibrators	101559	418305
Control 1	101560	418306
Control 2	101561	418306
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

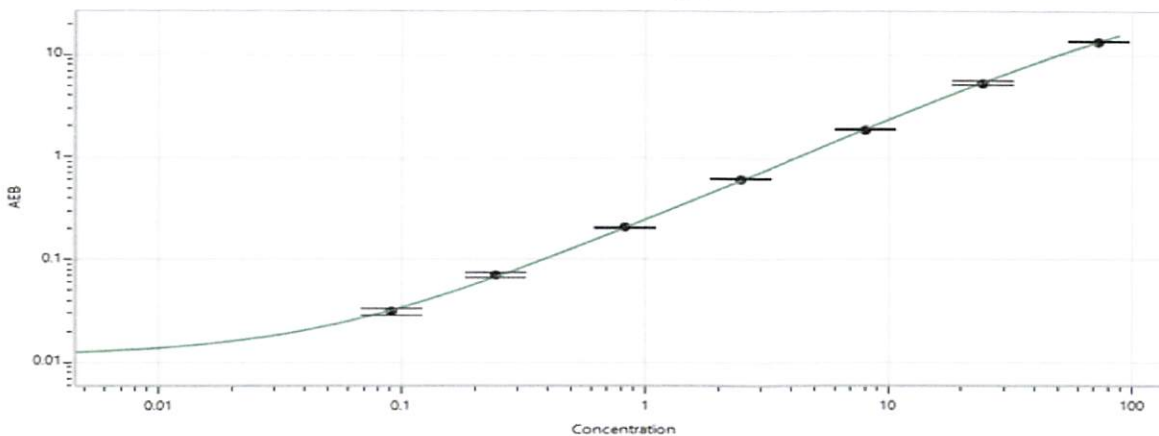
¹ Reagents are not Kit Lot Specific

† 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	418306	1.49	1.21 - 1.82
Control 2	418306	49.5	41.2 - 61.9

* 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	0.091
C	0.245
D	0.828
E	2.49
F	8.03
G	24.3
H	72.9

Review/Approval

Cody Carlson

QC Manager

[Signature] 07AUG2024

Name

Title

Signature/Date