

Product Number:	101552
Lot Number:	504721
Expiration Date:	15-Nov-26
Platform(s):	SR-X and HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	101554	513501
Detector Reagent	101553	525117
SBG Reagent	101555	513506
Sample Diluent Reagent	101556	513505
Reference Calibrators	101559	513210
Control 1	101560	513211
Control 2	101561	513211
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.

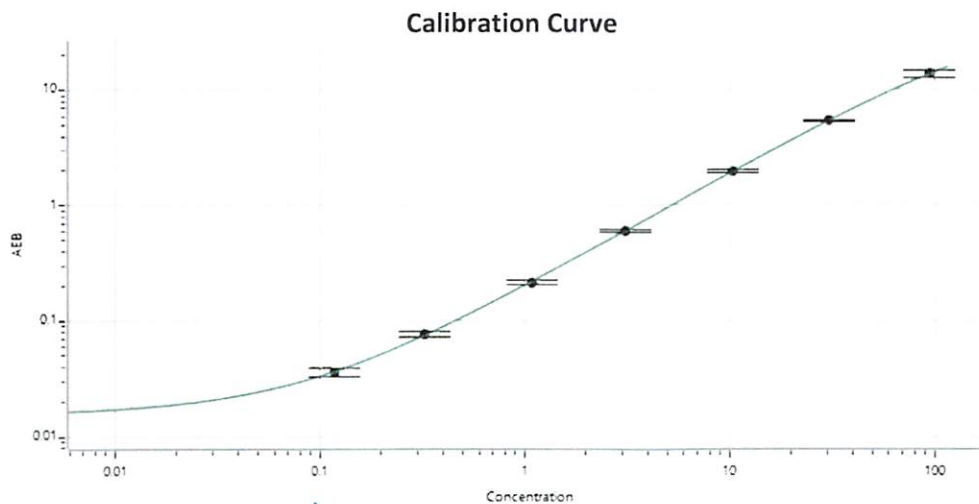
Refer to the Tau Advantage SR-X Kit Instructions (KI-0034).



Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	513211	0.959	0.864 - 1.30
Control 2	513211	59.2	47.3 - 74.7

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.



Calibrator Levels (pg/mL)

A	0.000
B	0.118
C	0.325
D	1.08
E	3.08
F	10.3
G	30.5
H	94.1

Review/Approval

Dhruvi Patel 8A Specialist II, 8A *SDP* 15 Sep 2025

Name

Title

Signature/Date