

Product Number:	101552
Lot Number:	504508
Expiration Date:	11-Aug-26
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	Control Lot Number
Bead Reagent	101554	507114	N/A
Detector Reagent	101553	509806	N/A
SBG Reagent	101555	504310	504310A
Sample Diluent Reagent	101556	504209	504209A
Reference Calibrators	101559	504504	504504A
Control 1	101560	504505	504505A
Control 2	101561	504505	504505A
SBG Diluent (SR-X)†	103158	N/A ¹	N/A ¹
RGP Reagent	103159	N/A ¹	N/A ¹

¹ Reagents are not Kit Lot Specific

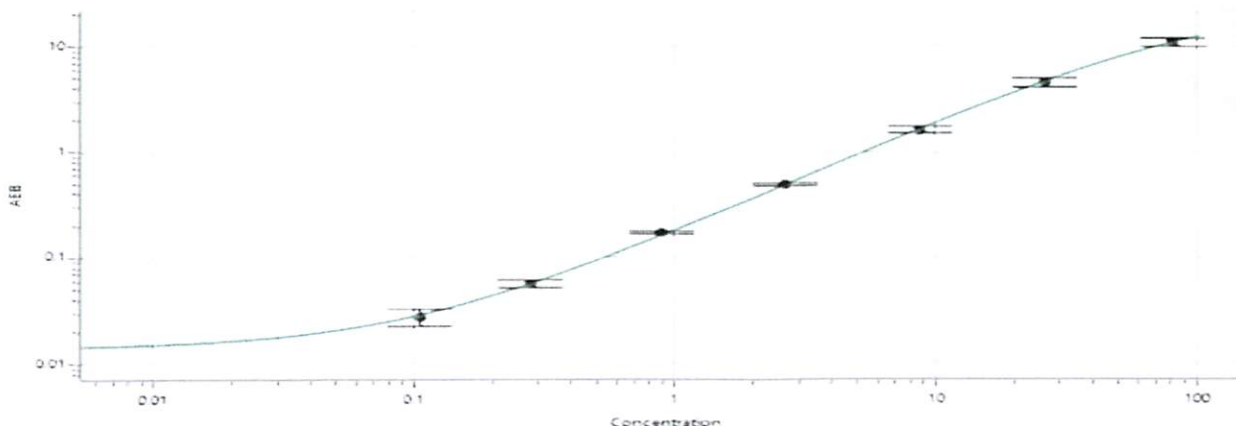
† Specific to use on the SR-X platform.

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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	504505	1.04	0.902 - 1.35
Control 2	504505	49.9	46.3 - 69.5

* 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	0.105
C	0.279
D	0.891
E	2.65
F	8.63
G	26.2
H	80.2

Review/Approval

Dipesh Patel Sr. Documental - QA *AM* 17 APR 2025

Name

Title

Signature/Date