

Product Number:	101552
Lot Number:	504720
Expiration Date:	17-Jan-27
Platform(s):	SR-X and HD-1/HD-X



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

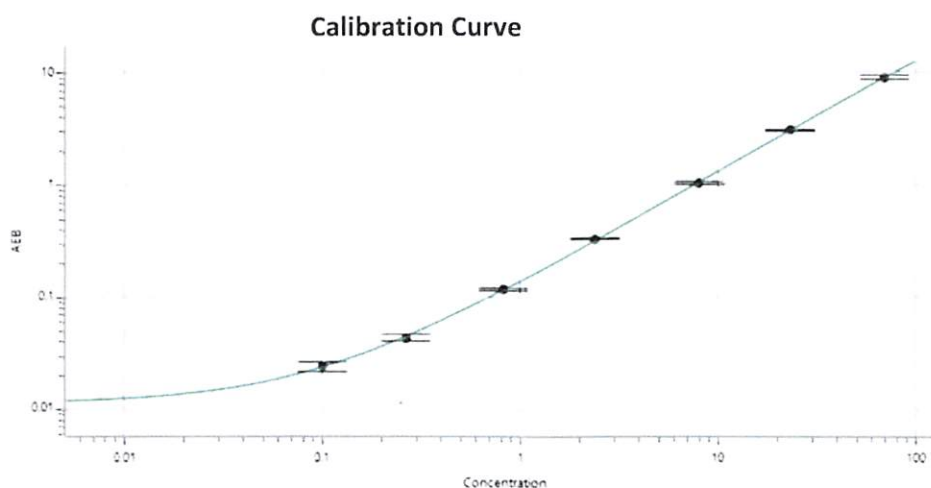
Component	Part Number	Lot Number
Bead Reagent	101554	519811
Detector Reagent	101553	525117
SBG Reagent	101555	519902
Sample Diluent Reagent	101556	519813
Reference Calibrators	101559	519701
Control 1	101560	519702
Control 2	101561	519702
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

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*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Spec. Range (pg/mL)	
Control 1	519702	0.885	0.785 - 1.37	
Control 2	519702	43.0	34.3 - 62.7	



Calibrator Levels (pg/mL)

A	0.000
B	0.100
C	0.264
D	0.819
E	2.37
F	7.93
G	23.0
H	68.9

Review/Approval

Dhruvi Patel QA Specialist II, QA DP/25SEP2025

Name

Title

Signature/Date

Quanterix Corporation

900 Middlesex Turnpike, Building 1, Billerica, 01821

techsupport@quanterix.com

©2025 Quanterix Corp. Simoa® is a registered trademark of Quanterix Corp.

For research use only. Not for use in diagnostic procedures.

Release

COA-0055 02