

Product Number:	101599
Lot Number:	504692
Expiration:	28-Oct-2026
Platform(s):	SR-X and HD-1/HD-X

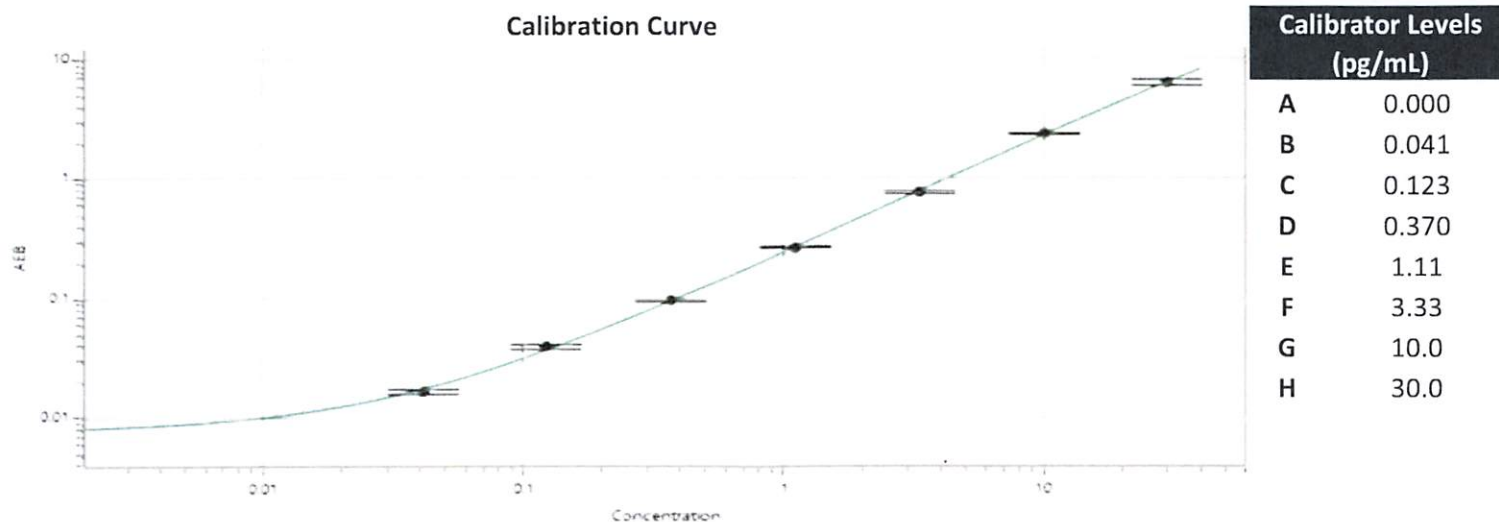
Component	Part Number	Lot Number
Bead Reagent	101601	522503
Detector Reagent	102012	525504
SBG Reagent	101602	522502
Sample Diluent	103886	522505
Calibrator Diluent	102014	511809
Calibrator Concentrate	100778	523308
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

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

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	523308	25.3	18.0 - 30.0	
Control 2	523308	0.561	0.375 - 0.625	

Calibrator Concentrate Value Assignment (pg/mL)	585
--	-----



Review/Approval

Dhruvi Patel QA Specialist II, QA GP / 22 Sep 2025

Name Title Signature/Date