

Product Number:	102039
Lot Number:	504708
Expiration:	28-Oct-2026
Platform(s):	SR-X and HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	102028	524602
Detector Stock	102029	523405
Calibrator Concentrate	102032	523807
Calibrator Diluent	102030	511810
Sample Diluent	103402	523409

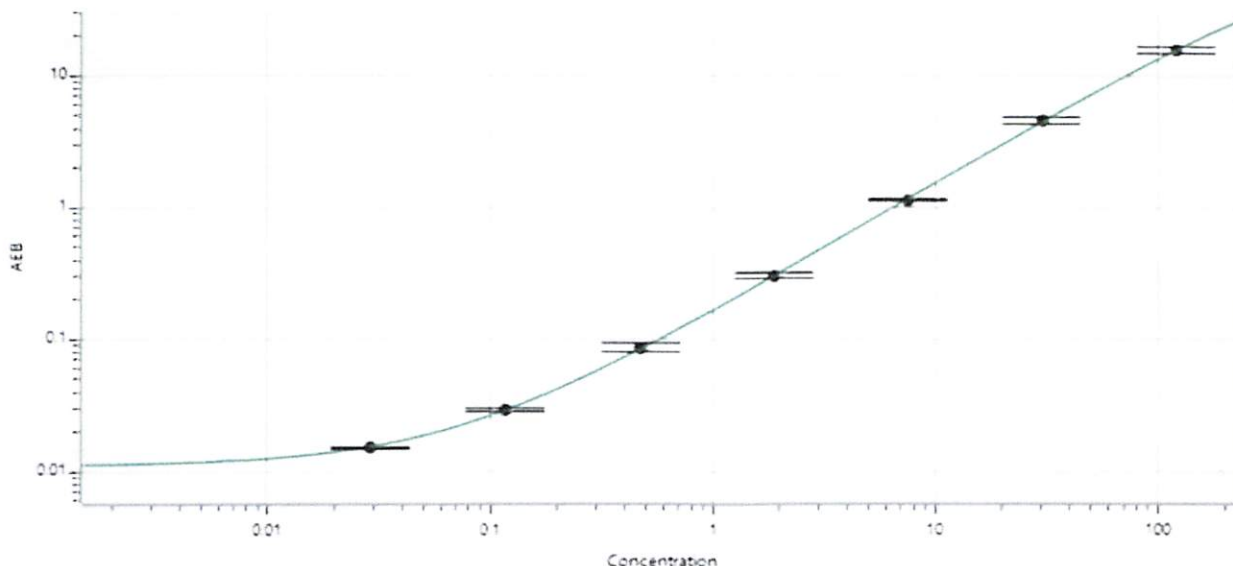
Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159



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

Calibrator Concentrate Target Value (ng/mL)	2.00
--	------

Calibration Curve



Review/Approval

Dhruvi Patel *QA Specialist II, QA* *DP* | 17 Sep 2025

Name

Title

Signature/Date