

Product Number:	102215
Lot Number:	536639
Expiration:	29-Mar-2027
Platform(s):	SR-X and HD-1/HD-X

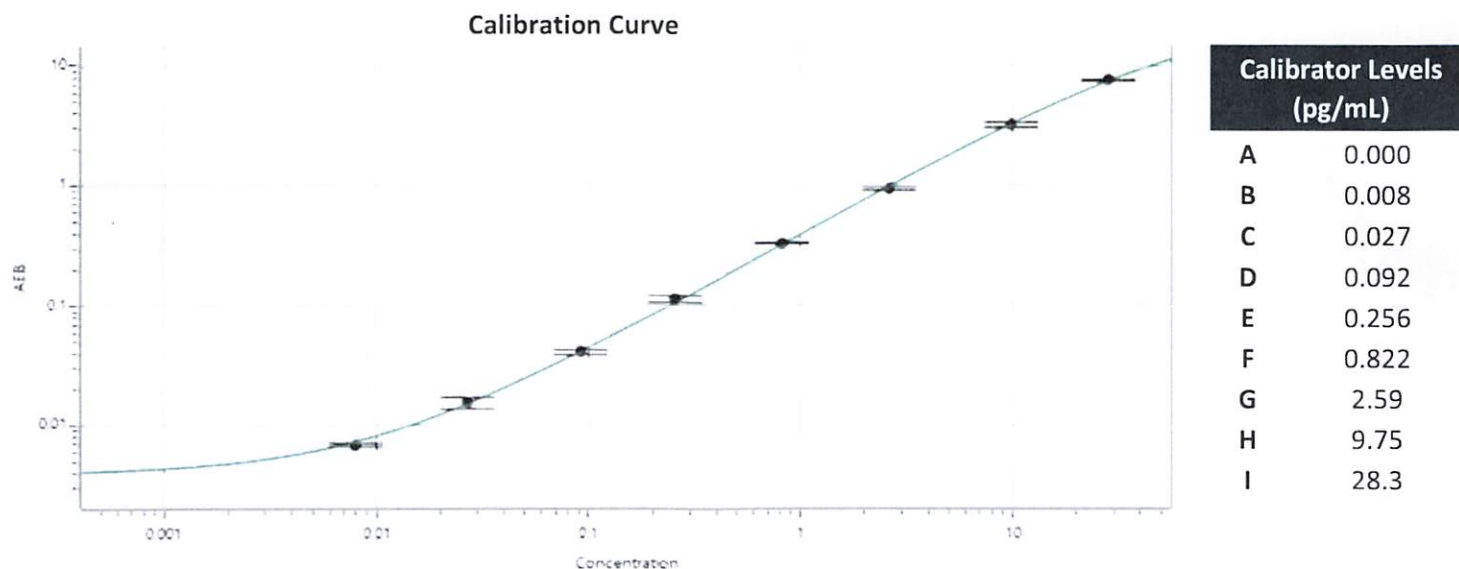
Component	Part Number	Lot Number
Bead Reagent	104297	531003
Detector Reagent	104338	531007
SBG Reagent	104340	531005
Sample Diluent	104342	531008
Calibrators	104343	531405
Control 1	104344	531406
Control 2	104345	531406
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	531406	0.088	0.071 - 0.107
Control 2	531406	9.15	7.26 - 11.4

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



Review/Approval

Dhruvi Patel QA Specialist II DP / 04 Dec 2025

Name Title Signature/Date