

Product Number:	105569
Lot Number:	504447
Expiration:	4-May-2026
Platform(s):	HD-X



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

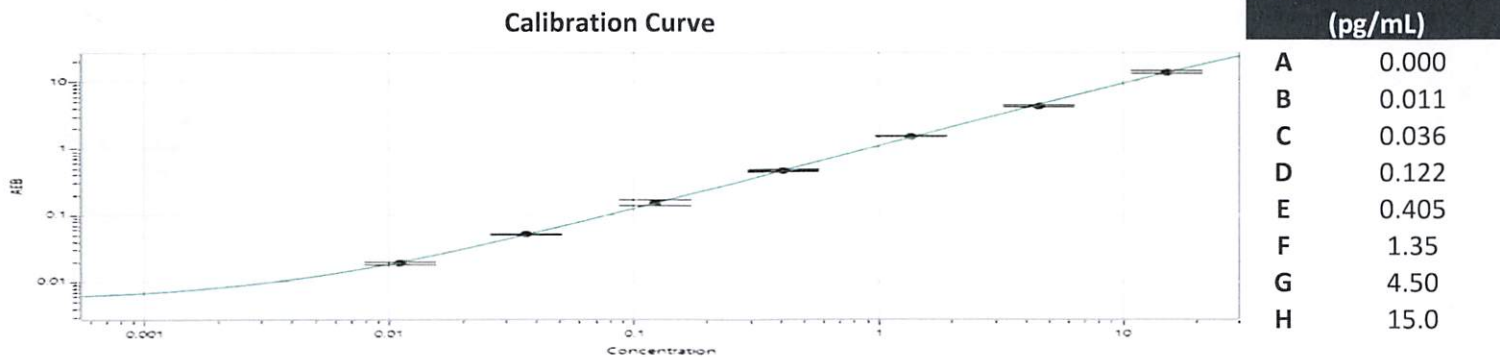
Component	Part Number	Lot Number
Bead Reagent	105548	432503
Detector Reagent	105550	435301
SBG Reagent	105552	432602
Sample Diluent	105553	430924
Calibrator Diluent	105555	430926
Control Diluent	105556	430925
Calibrator Concentrate	105565	427521
Control 1 Concentrate	105567	427523
Control 2 Concentrate	105568	427524
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	427523	0.708	0.529 - 0.983
Control 2 Concentrate	427524	10.4	8.30 - 15.4

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

Calibrator Concentrate Reconstitution Volume (uL)	1072
Control 1 Concentrate Reconstitution Volume (uL)	1072
Control 2 Concentrate Reconstitution Volume (uL)	1072



Review/Approval Dhruvi Patel DOC Control I, QA SDP 28 Jan 2025
 Name Title Signature/Date