

Product Number:	106068
Lot Number:	504679
Expiration:	2-Jan-2027
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Bead Reagent	104546	520403
Detector Reagent	104548	520504
SBG Reagent	104550	520505
Sample Diluent	104552	518311
Reference Calibrators	106071	519507
Control 1	106069	519508
Control 2	106070	519509
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

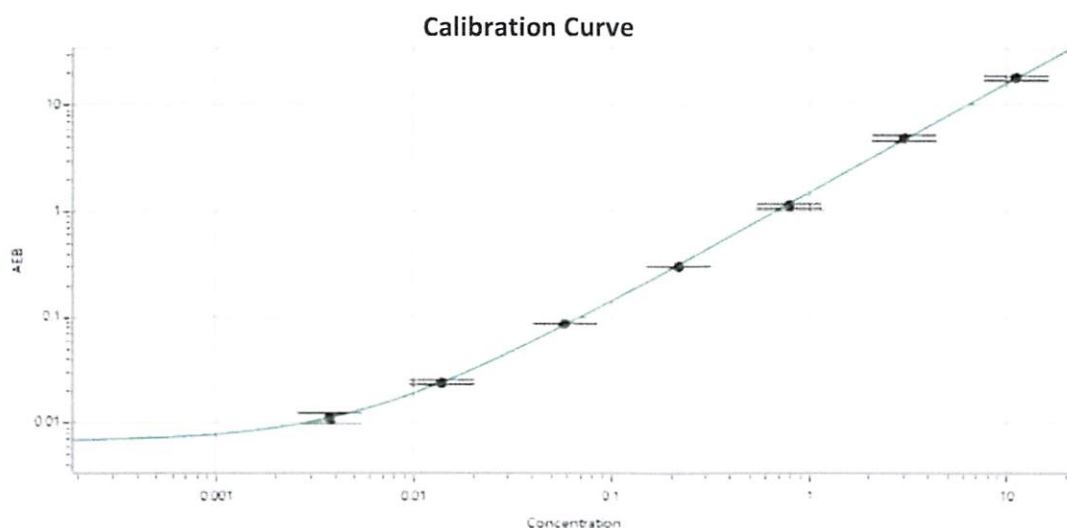
¹ Reagents are not Kit Lot Specific



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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	519508	0.665	0.445 - 0.826
Control 2	519509	6.69	4.55 - 8.45

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



Review/Approval

Dhruvi Patel QA Specialist II, QA SDP / 15 Sep 2025

Name

Title

Signature/Date