

Product Number:	105827
Lot Number:	536600
Expiration:	17-Jan-2027
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	105803	521306
Detector Reagent	105805	527505
SBG Reagent	105807	521705
Sample Diluent	105808	519818
Calibrator Diluent	105810	519820
Control Diluent	105811	519819
Calibrator Concentrate	105823	504914
Control 1 Concentrate	105825	504916
Control 2 Concentrate	105826	504917
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

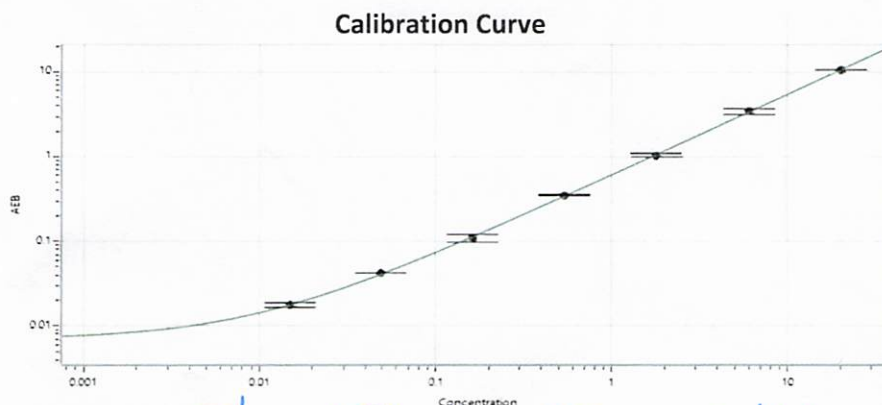


Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	504916	0.427	0.329 - 0.611
Control 2 Concentrate	504917	9.76	6.99 - 13.0

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

Calibrator Concentrate Reconstitution Volume (uL)	919
Control 1 Concentrate Reconstitution Volume (uL)	919
Control 2 Concentrate Reconstitution Volume (uL)	919



Calibrator Levels (pg/mL)	
A	0.000
B	0.015
C	0.049
D	0.162
E	0.540
F	1.80
G	6.00
H	20.0

Review/Approval

Dhwani Patel
Name

QA Specialist II
Title

SDP / 17 Nov 2025
Signature/Date