

Product Number:	105634
Lot Number:	504709
Expiration:	27-Feb-2027
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	105608	524601
Detector Reagent	105610	524710
SBG Reagent	105612	524712
Sample Diluent	105614	523910
Calibrator Diluent	105617	523912
Control Diluent	105618	523911
Calibrator Concentrate	105630	430514
Control 1 Concentrate	105632	430516
Control 2 Concentrate	105633	430517
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific



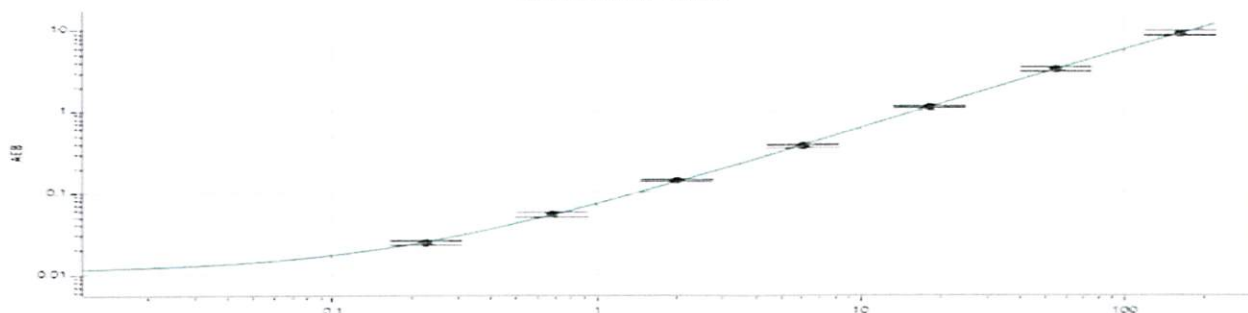
Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	430516	7.56	5.70 - 10.6
Control 2 Concentrate	430517	116	88.4 - 164

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

Calibrator Concentrate Reconstitution Volume (uL)	746
Control 1 Concentrate Reconstitution Volume (uL)	746
Control 2 Concentrate Reconstitution Volume (uL)	746

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.225
C	0.675
D	2.02
E	6.07
F	18.2
G	54.7
H	164

Dipesh Patel Sr. QA Specialist [Signature] 26 Feb 2027
 Name Title Signature/Date