

Product Number:	105187
Lot Number:	536628
Expiration:	17-Jan-2027
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	105139	529604
Detector Reagent	105141	530107
SBG Reagent	105143	530108
Sample Diluent	105144	519821
Calibrator Diluent	105172	528828
Control Diluent	105173	528829
Calibrator Concentrate	105183	412323
Control 1 Concentrate	105185	412324
Control 2 Concentrate	105186	412325
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific



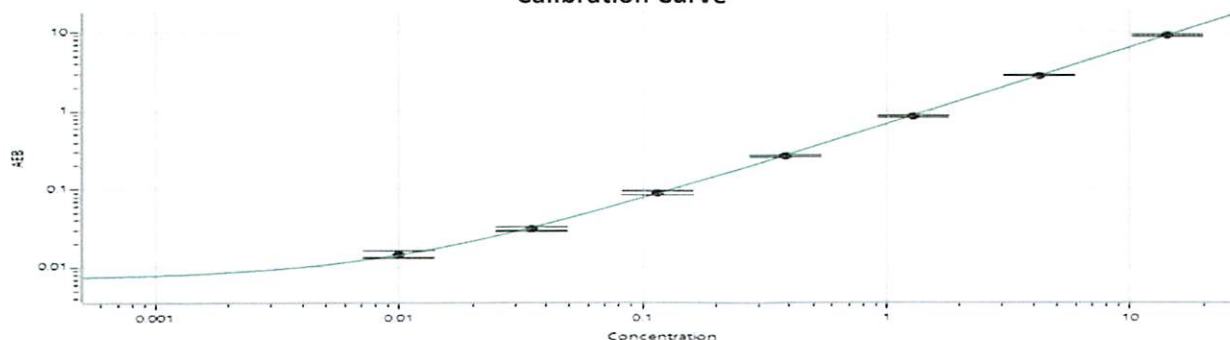
Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412324	0.726	0.517 - 0.960
Control 2 Concentrate	412325	17.2	12.4 - 22.9

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

Calibrator Concentrate Reconstitution Volume (uL)	918
Control 1 Concentrate Reconstitution Volume (uL)	918
Control 2 Concentrate Reconstitution Volume (uL)	918

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.010
C	0.035
D	0.115
E	0.385
F	1.28
G	4.26
H	14.2

Review/Approval

Dhruvi Patel *QA Specialist II* *DP/09Dec2025*
 Name Title Signature/Date