

Product Number:	104364
Lot Number:	504688
Expiration:	5-Oct-2027
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Bead Reagent	104352	523006
Detector Reagent	104353	523005
SBG Reagent	104354	521703
Sample Diluent Reagent	104355	521801
Calibrator Diluent Reagent	104358	521716
Control Diluent Reagent	104359	521717
Calibrator Concentrate	104361	412713
Control 1 Concentrate	104362	412714
Control 2 Concentrate	104363	412715
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific



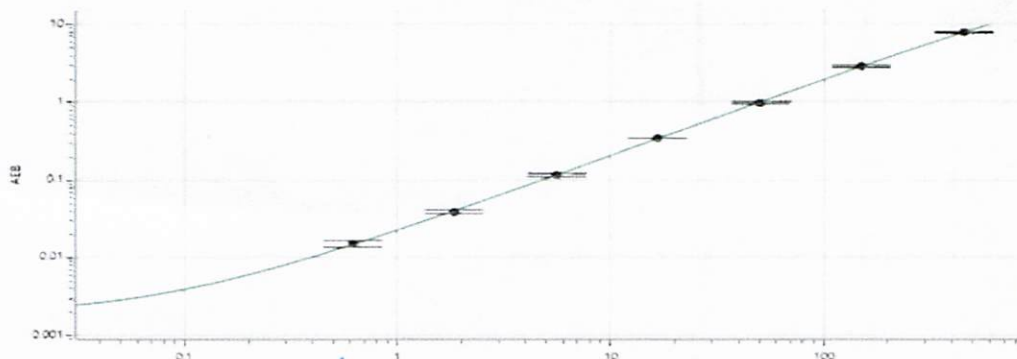
Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	412714	4.43	2.84 - 7.52
Control 2	412715	324	282 - 562

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

Calibrator Concentrate Reconstitution Volume (uL)	1500
Control 1 Concentrate Reconstitution Volume (uL)	1500
Control 2 Concentrate Reconstitution Volume (uL)	1500

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.617
C	1.85
D	5.56
E	16.7
F	50.0
G	150
H	450

Review/Approval

Dhwani Patel

Name

QA Specialist II

Title

DP / 26 Nov 2025

Signature/Date