

Product Number:	104428
Lot Number:	504498
Expiration:	19-Feb-2026
Platform(s):	HD-X and SR-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	104409	503703	503703A
Detector Reagent	104411	504202	504202A
SBG Reagent	104413	504203	504203A
Sample Diluent	104415	503510	503510A
Calibrator Diluent	104418	423225	N/A
Control Diluent	104476	423226	N/A
Calibrator Concentrate	104483	412912	N/A
Control 1 Concentrate	104426	412910	N/A
Control 2 Concentrate	104427	412911	N/A
RGP 2.0 Reagent	104321	N/A ¹	N/A ¹
Activation Buffer Reagent	104324	N/A ¹	N/A ¹

¹ Reagents are not Kit Lot Specific

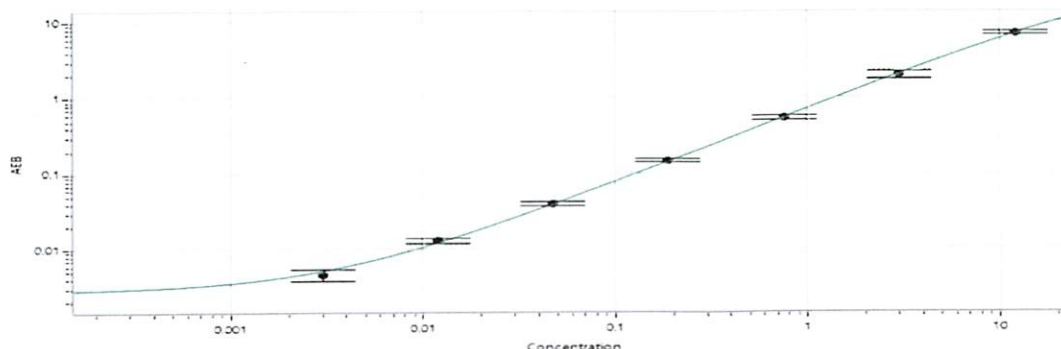
Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Spec. Range (pg/mL)	
Control 1 Concentrate	412910	0.188	0.099 - 0.241	
Control 2 Concentrate	412911	5.28	3.76 - 6.98	

Calibrator Concentrate Reconstitution Volume (uL)	893
Control 1 Concentrate Reconstitution Volume (uL)	893
Control 2 Concentrate Reconstitution Volume (uL)	893

Calibrator Levels (pg/mL)

A	0.000
B	0.003
C	0.012
D	0.047
E	0.188
F	0.750
G	3.00
H	12.0

Calibration Curve



Review/Approval

Dipesh Patel Sr. Doc control
Name Title

[Signature] 13 MAR 2025
Signature/Date