

Product Number:	104542
Lot Number:	504497
Expiration:	10-Aug-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	104515	509409	N/A
Detector Reagent	104517	504301	504301A
SBG Reagent	104519	504302	N/A
Sample Diluent	104521	504124	504124A
Calibrator Diluent	104524	506317	N/A
Control Diluent	104525	506318	N/A
Calibrator Concentrate	104538	400911	N/A
Control 1 Concentrate	104540	400912	N/A
Control 2 Concentrate	104541	400913	N/A
RGP 2.0 Reagent	104321	N/A ¹	N/A ¹
Activation Buffer Reagent	104324	N/A ¹	N/A ¹

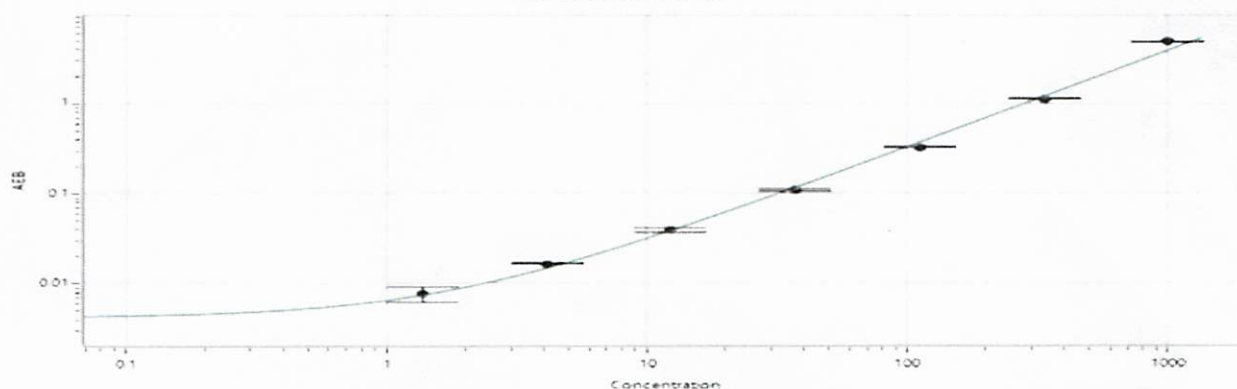
¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	400912	377	264 - 491
Control 2 Concentrate	400913	4484	3201 - 5944

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

Calibrator Concentrate Reconstitution Volume (uL)	1177
Control 1 Concentrate Reconstitution Volume (uL)	975
Control 2 Concentrate Reconstitution Volume (uL)	975

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Dilip Patel Sr. Doc Control - QA TM / 25 Jun 2025
Name Title Signature/Date