

Product Number:	105737
Lot Number:	504577
Expiration:	27-Sep-2026
Platform(s):	HD-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	105712	510007	N/A
Detector Reagent	105714	512504	N/A
SBG Reagent	105716	512505	N/A
Sample Diluent	105718	508609	N/A
Calibrator Diluent	105720	511506	N/A
Control Diluent	105721	511508	N/A
Calibrator Concentrate	105733	504410	504410A
Control 1 Concentrate	105735	504412	504412A
Control 2 Concentrate	105736	504413	504413A
RGP 2.0 Reagent	104321	N/A ¹	N/A ¹
Activation Buffer Reagent	104324	N/A ¹	N/A ¹

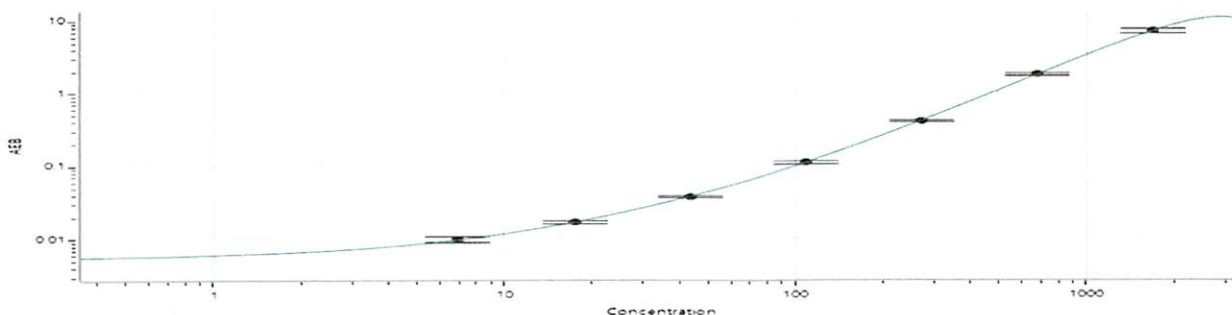
¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	504412A	2066	1647 - 3059
Control 2 Concentrate	504413A	20798	17132 - 31817

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

Calibrator Concentrate Reconstitution Volume (uL)	762
Control 1 Concentrate Reconstitution Volume (uL)	762
Control 2 Concentrate Reconstitution Volume (uL)	300

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	6.96
C	17.4
D	43.5
E	109
F	272
G	680
H	1700

Review/Approval

Dhruvi Patel Doc Control II, QA *DP* / 30 Jun 2025

Name

Title

Signature/Date