

Product Number:	104570
Lot Number:	504614
Expiration:	21-Sep-2026
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Bead Reagent	104546	508010
Detector Reagent	104548	508401
SBG Reagent	104550	508402
Sample Diluent	104552	508417
Calibrator Diluent Reagent	104745	508416
Control Diluent Reagent	104554	508415
Calibrator Concentrate	104566	405210
Control 1 Concentrate	104568	405211
Control 2 Concentrate	104569	405212
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.

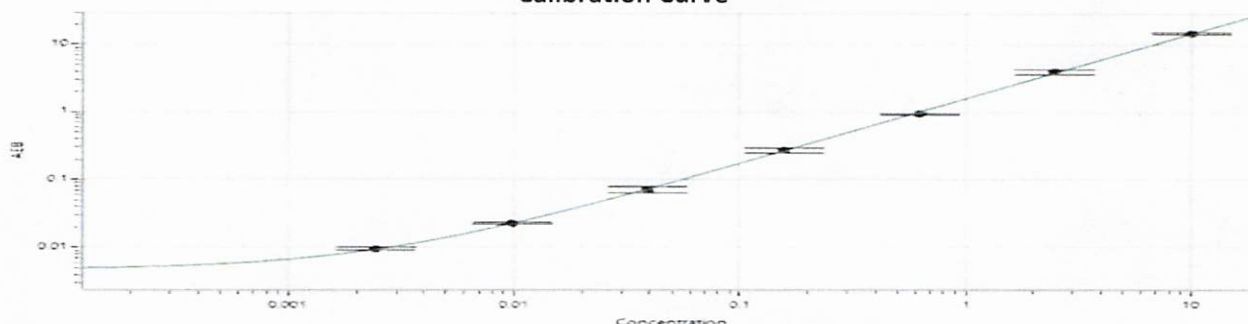
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	405211	0.748	0.516 - 0.958
Control 2 Concentrate	405212	7.64	5.40 - 10.0

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

Calibrator Concentrate Reconstitution Volume (uL)	976
Control 1 Concentrate Reconstitution Volume (uL)	976
Control 2 Concentrate Reconstitution Volume (uL)	976

Calibrator Levels (pg/mL)	
A	0.000
B	0.00244
C	0.00977
D	0.0391
E	0.156
F	0.625
G	2.50
H	10.0

Calibration Curve



Review/Approval

Dipesh Patel

Name

Sr. Doc. Control - QA

Title

ET / 02 Jun 2025

Signature/Date