

Product Number:	104570
Lot Number:	504525
Expiration:	10-Sep-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	104546	507013	N/A
Detector Reagent	104548	507108	N/A
SBG Reagent	104550	507109	N/A
Sample Diluent	104552	506915	N/A
Calibrator Diluent Reagent	104745	506914	506914A
Control Diluent Reagent	104554	506913	N/A
Calibrator Concentrate	104566	405210	N/A
Control 1 Concentrate	104568	405211	N/A
Control 2 Concentrate	104569	405212	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

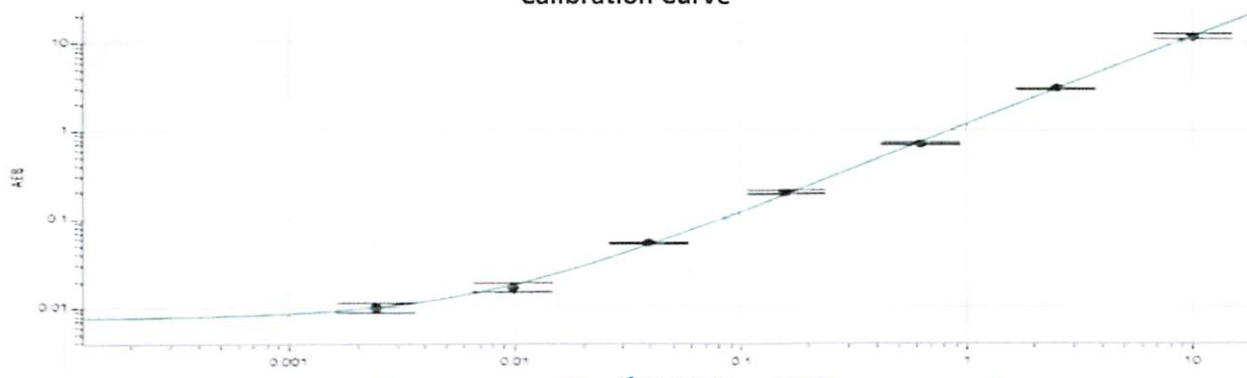
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	405211	0.758	0.532 - 0.987
Control 2 Concentrate	405212	8.28	4.94 - 10.2

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

Calibrator Concentrate Reconstitution Volume (uL)	1122
Control 1 Concentrate Reconstitution Volume (uL)	1122
Control 2 Concentrate Reconstitution Volume (uL)	1122

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 *Sr. Documental-RA* *em 1/16 APR 2025*
 Name Title Signature/Date