

Product Number:	104570
Lot Number:	504683
Expiration:	15-Feb-2027
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Bead Reagent	104546	523007
Detector Reagent	104548	524806
SBG Reagent	104550	524807
Sample Diluent	104552	522709
Calibrator Diluent Reagent	104745	522711
Control Diluent Reagent	104554	522710
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	0.726	0.503 - 0.934
Control 2 Concentrate	405212	6.93	5.00 - 9.29

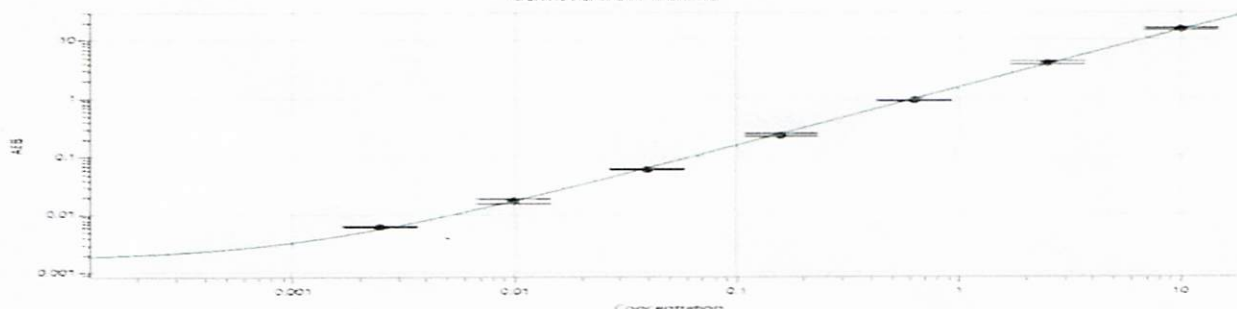
* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	877
Control 1 Concentrate Reconstitution Volume (uL)	877
Control 2 Concentrate Reconstitution Volume (uL)	877

Calibrator Levels (pg/mL)

A	0
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

Dilip Patel

Name

Sr. QASpecialist

Title

29 SEP 2025

Signature/Date