

Product Number:	104570
Lot Number:	504576
Expiration:	13-Nov-2026
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Bead Reagent	104546	513302
Detector Reagent	104548	513503
SBG Reagent	104550	513504
Sample Diluent	104552	513305
Calibrator Diluent Reagent	104745	513307
Control Diluent Reagent	104554	513306
Calibrator Concentrate	104566	406408
Control 1 Concentrate	104568	406409
Control 2 Concentrate	104569	406410
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.

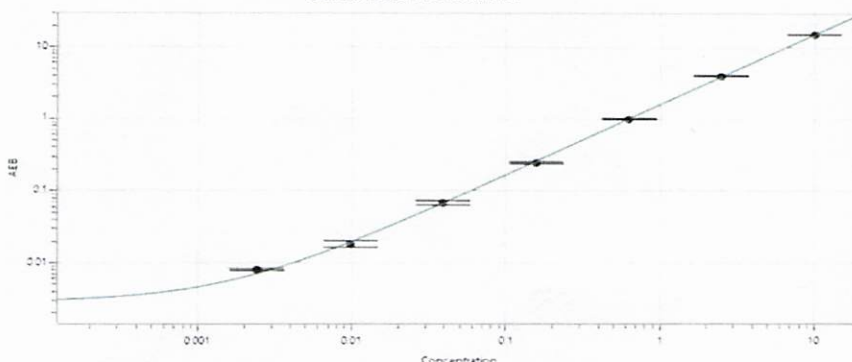
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	406409	0.736	0.537 - 0.998
Control 2 Concentrate	406410	8.36	5.27 - 9.79

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

Calibrator Concentrate Reconstitution Volume (uL)	868
Control 1 Concentrate Reconstitution Volume (uL)	868
Control 2 Concentrate Reconstitution Volume (uL)	868

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

Diipul Patel *Sr. Doc. control-QA* *ET* / 16 Jun 2025
 Name Title Signature/Date