

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

Component	Part Number	Lot Number
Bead Reagent	104546	516101
Detector Reagent	104548	516303
SBG Reagent	104550	516304
Sample Diluent	104552	516105
Calibrator Diluent Reagent	104745	516107
Control Diluent Reagent	104554	516106
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

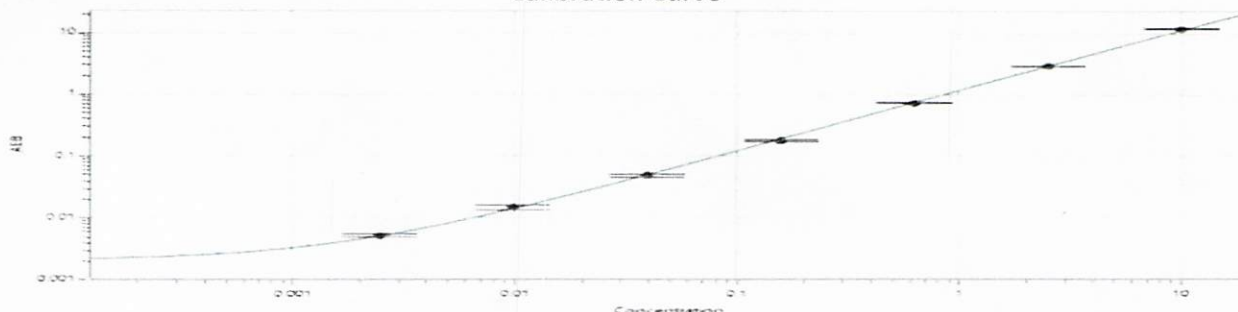
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	406409	0.672	0.520 - 0.966
Control 2 Concentrate	406410	7.51	5.40 - 10.0

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

Calibrator Concentrate Reconstitution Volume (uL)	963
Control 1 Concentrate Reconstitution Volume (uL)	963
Control 2 Concentrate Reconstitution Volume (uL)	963

**Calibrator Levels
(pg/mL)**

Calibration Curve



A	0
B	0.00244
C	0.00977
D	0.0391
E	0.156
F	0.625
G	2.50
H	10.0

Review/Approval

Dipesh Patel
Name

Ser. Doc. Control-QA

Signature/Date