

Product Number:	104570
Lot Number:	504482
Expiration:	6-Sep-2025
Platform(s):	HD-X and SR-X



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

Component	Part Number	Lot Number
Bead Reagent	104546	422503
Detector Reagent	104548	422712
SBG Reagent	104550	422713
Sample Diluent	104552	406619
Calibrator Diluent Reagent	104745	422621
Control Diluent Reagent	104554	422620
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

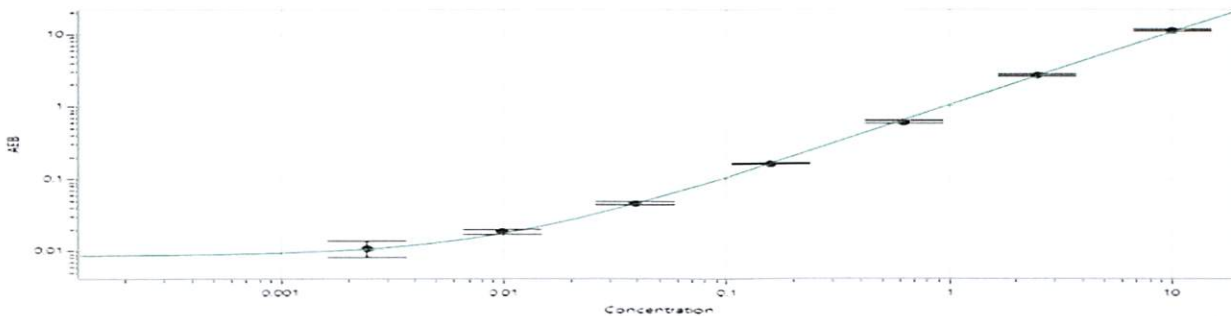
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	406409	0.823	0.449 - 0.834
Control 2 Concentrate	406410	6.98	4.41 - 8.19

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

Calibrator Concentrate Reconstitution Volume (uL)	972
Control 1 Concentrate Reconstitution Volume (uL)	972
Control 2 Concentrate Reconstitution Volume (uL)	972

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

Doc Control II RA
Title

et 18 DEC 2024
Signature/Date