

Product Number:	104364
Lot Number:	504501
Expiration:	7-Aug-2026
Platform(s):	HD-X and SR-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	104352	502807	503807A
Detector Reagent	104353	416212	N/A
SBG Reagent	104354	420509	N/A
Sample Diluent Reagent	104355	416201	N/A
Calibrator Diluent Reagent	104358	415913	N/A
Control Diluent Reagent	104359	415914	N/A
Calibrator Concentrate	104361	404519	N/A
Control 1 Concentrate	104362	404520	N/A
Control 2 Concentrate	104363	404521	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

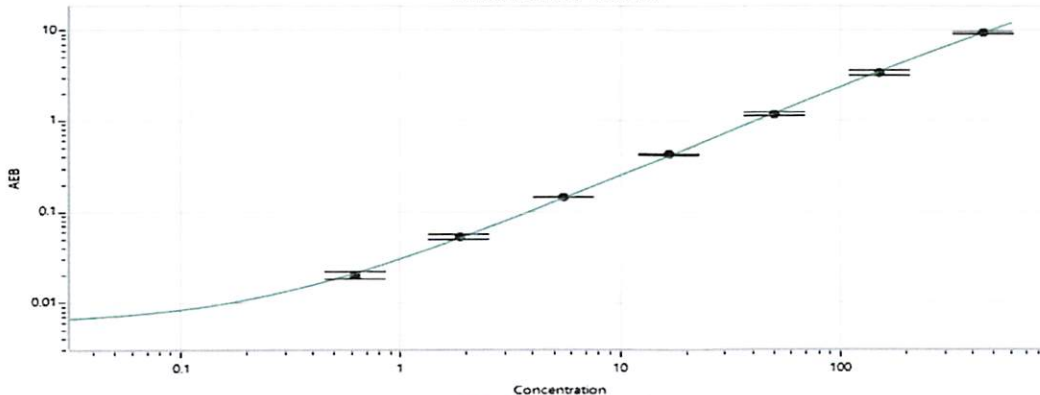
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	404520	3.85	3.32 - 6.25
Control 2	404521	324	272 - 506

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

Calibrator Concentrate Reconstitution Volume (uL)	1387
Control 1 Concentrate Reconstitution Volume (uL)	1387
Control 2 Concentrate Reconstitution Volume (uL)	1387

Calibrator Levels (pg/mL)	
A	0
B	0.617
C	1.85
D	5.56
E	16.7
F	50.0
G	150
H	450

Calibration Curve



Review/Approval: Dipesh Patel / Doc control TQA / [Signature] / 16 Feb 2025

Name: Dipesh Patel Title: Doc control TQA Signature/Date: [Signature] / 16 Feb 2025