

Product Number:	104364
Lot Number:	504506
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	503101	503101A
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

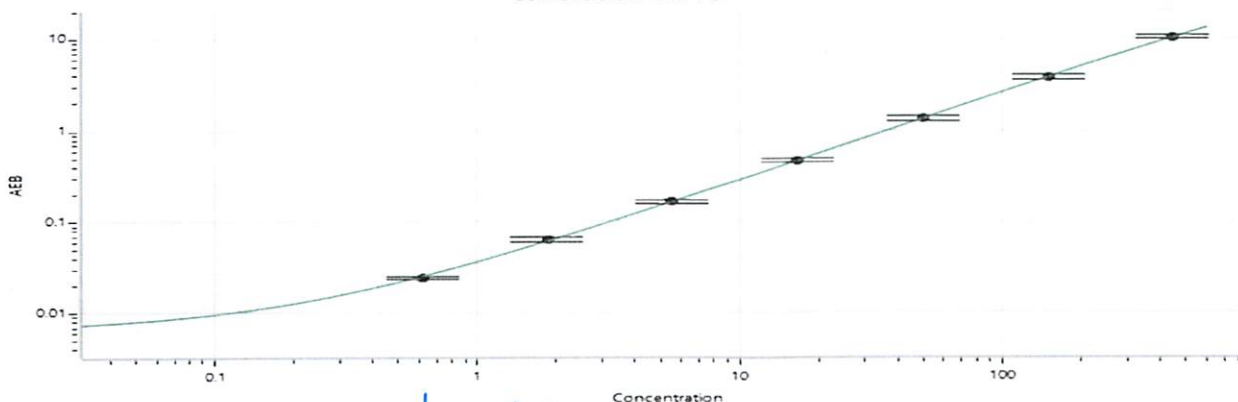
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	404520	3.66	2.86 - 6.37
Control 2	404521	332	278 - 516

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

Calibrator Concentrate Reconstitution Volume (uL)	1401
Control 1 Concentrate Reconstitution Volume (uL)	1401
Control 2 Concentrate Reconstitution Volume (uL)	1401

Calibrator Levels (pg/mL)	
A	0.000
B	0.617
C	1.85
D	5.56
E	16.7
F	50.0
G	150
H	450

Calibration Curve



Review/Approval

Dhwani Patel *DOC Control II, QA* *SDP* *04 Apr 2025*

Name

Title

Signature/Date