

Product Number:	104618
Lot Number:	504238
Expiration:	23-Feb-2026
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	104602	414402
Detector Reagent	104603	415003
SBG Reagent	104604	423605
Sample Diluent	104605	414403
Calibrator Diluent	104607	414414
Control Diluent	104608	414415
Calibrator Concentrate	104614	400907
Control 1 Concentrate	104616	400908
Control 2 Concentrate	104617	400909
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

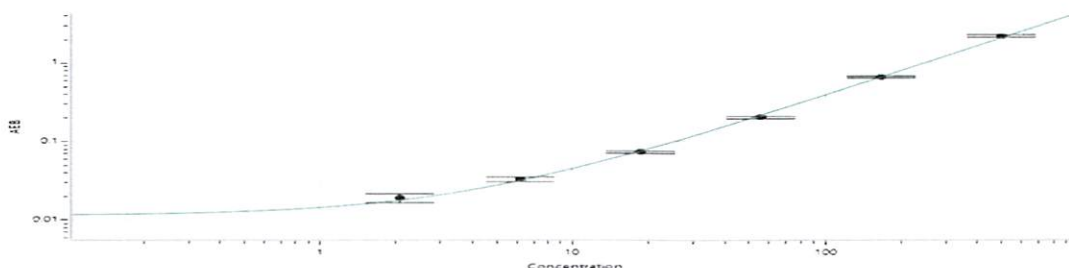
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	400908	66.0	43.8 - 85.1
Control 2 Concentrate	400909	1865	1167 - 2167

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

Calibrator Concentrate Reconstitution Volume (uL)	1517
Control 1 Concentrate Reconstitution Volume (uL)	841
Control 2 Concentrate Reconstitution Volume (uL)	841

Calibrator Levels (pg/mL)	
A	0.000
B	2.06
C	6.17
D	18.5
E	55.6
F	167
G	500

Calibration Curve



Review/Approval

TAM LEE DOC CONTROL II, QA Tam Lee 15 Nov 2024

Name

Title

Signature/Date