

Product Number:	105605
Lot Number:	504449
Expiration:	4-May-2026
Platform(s):	HD-X



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

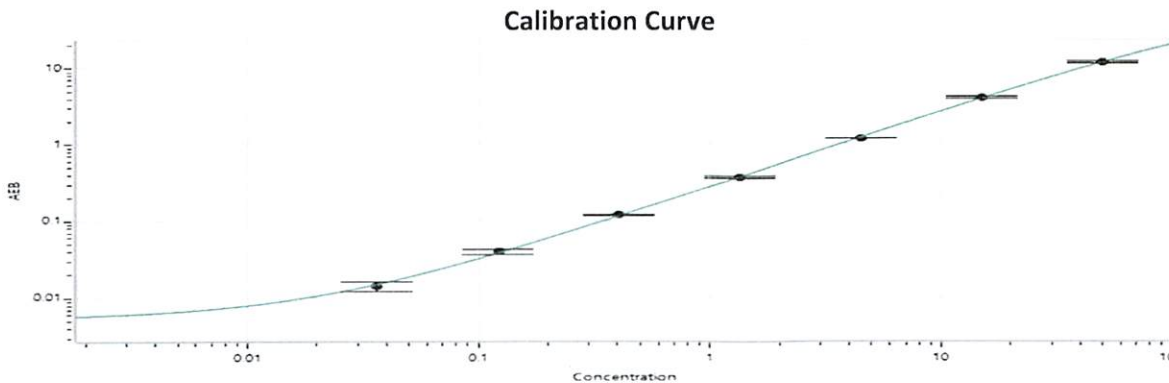
Component	Part Number	Lot Number
Bead Reagent	105593	432707
Detector Reagent	105595	433701
SBG Reagent	105597	433702
Sample Diluent	105599	432005
Calibrator Diluent	105601	430922
Control Diluent	105602	430923
Calibrator Concentrate	105522	427513
Control 1 Concentrate	105524	427515
Control 2 Concentrate	105525	427516
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	427515	1.35	0.801 - 1.95
Control 2 Concentrate	427516	85.8	68.9 - 128

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

Calibrator Concentrate Reconstitution Volume (uL)	1183
Control 1 Concentrate Reconstitution Volume (uL)	1183
Control 2 Concentrate Reconstitution Volume (uL)	1183



Calibrator Levels (pg/mL)	
A	0.000
B	0.036
C	0.122
D	0.405
E	1.35
F	4.50
G	15.0
H	50.0

Review/Approval Dhwani Patel DOC Control I, QA DP 31 Dec 2024
 Name Title Signature/Date