

Product Number:	104512
Lot Number:	504321
Expiration:	23-Jan-2026
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	104485	420505
Detector Reagent	104487	421306
SBG Reagent	104489	421307
Sample Diluent	104491	420510
Calibrator Diluent	104494	420512
Control Diluent	104495	420511
Calibrator Concentrate	104508	330613
Control 1 Concentrate	104510	330615
Control 2 Concentrate	104511	330616
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

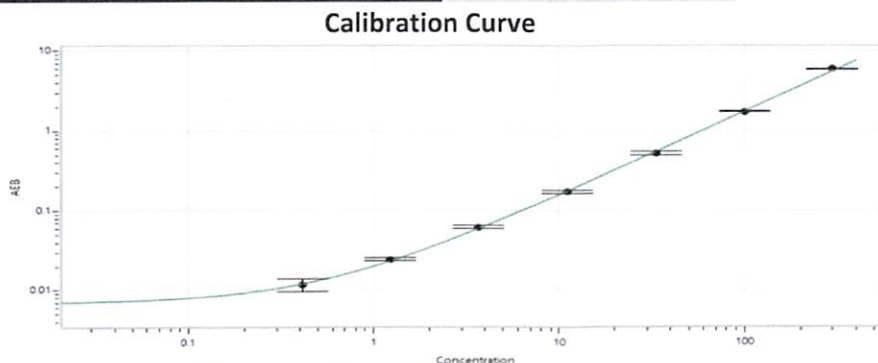
¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	330615	3.46	2.41 - 4.48
Control 2 Concentrate	330616	265	175 - 325

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

Calibrator Concentrate Reconstitution Volume (uL)	811
Control 1 Concentrate Reconstitution Volume (uL)	811
Control 2 Concentrate Reconstitution Volume (uL)	811

Calibrator Levels (pg/mL)	
A	0
B	0.412
C	1.23
D	3.70
E	11.1
F	33.3
G	100
H	300



Review/Approval

Dhvani Patel
Name

DOC Control I QA
Title

DPP / 12 Sep 2024
Signature/Date