

Product Number:	104588
Lot Number:	504407
Expiration:	12-Jan-2026
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	104572	427105
Detector Reagent	104573	427401
SBG Reagent	104574	427402
Sample Diluent Reagent	104575	426211
Calibrator Diluent Reagent	104577	419407
Control Diluent Reagent	104578	419408
Calibrator Concentrate	104584	329610
Control 1 Concentrate	104586	329612
Control 2 Concentrate	104587	329613
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

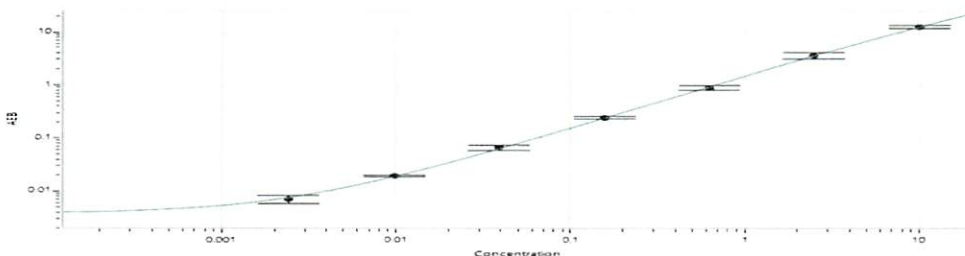
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	329612	0.103	0.068 - 0.126
Control 2 Concentrate	329613	2.47	1.75 - 3.25

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

Calibrator Concentrate Reconstitution Volume (uL)	806
Control 1 Concentrate Reconstitution Volume (uL)	806
Control 2 Concentrate Reconstitution Volume (uL)	806

Calibrator Levels (pg/mL)	
---------------------------	--

Calibration Curve



A	0
B	0.002
C	0.010
D	0.039
E	0.156
F	0.625
G	2.50
H	10.0

Review/Approval

Dhruvi Patel *DOC Control, I, QA* *DP* *08 Nov 2024*
Name Title Signature/Date