

Component	Part Number	Lot Number
Calibrator Concentrate	104466	317333
Control 1 Concentrate	104468	402941
Control 2 Concentrate	104469	402942



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

Calibrator Concentrate Reconstitution Volume (uL)	1144
Control 1 Concentrate Reconstitution Volume (uL)	1144
Control 2 Concentrate Reconstitution Volume (uL)	1144

Review/Approval

Dhruvi Patel QA Specialist II, QA DP / 26 Aug 2025

Name Title Signature/Date