

Component	Part Number	Lot Number
Calibrator Concentrate	104566	405210
Control 1 Concentrate	104568	405211
Control 2 Concentrate	104569	405212



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

Calibrator Concentrate Reconstitution Volume (uL)	877
Control 1 Concentrate Reconstitution Volume (uL)	877
Control 2 Concentrate Reconstitution Volume (uL)	877

Review/Approval

Dhruvi Patel QA Specialist II [Signature] / 26 Nov 2025
Name Title Signature/Date