

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

Calibrator Concentrate Reconstitution Volume (uL)	1393
Control 1 Concentrate Reconstitution Volume (uL)	1393
Control 2 Concentrate Reconstitution Volume (uL)	1393

Review/Approval

Dhvani Patel QA Specialist II DP / 04 Dec 2025

Name Title Signature/Date