

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

Calibrator Concentrate Reconstitution Volume (uL)	1250
Control 1 Concentrate Reconstitution Volume (uL)	1250
Control 2 Concentrate Reconstitution Volume (uL)	1250

William Chan	document control Specialist I	William Chan 04DEC2025
Name	Title	Signature/Date