

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

Calibrator Concentrate Reconstitution Volume (uL)	972
Control 1 Concentrate Reconstitution Volume (uL)	972
Control 2 Concentrate Reconstitution Volume (uL)	972

Review/Approval

Dhvani Patel

DOC Control II, QA

GP 01 May 2025

Name

Title

Signature/Date

Quanterix Corporation
900 Middlesex Turnpike, Building 1, Billerica, 01821
techsupport@quanterix.com

©2025 Quanterix Corp. Simoa® is a registered trademark of Quanterix Corp.
1 For research use only. Not for use in diagnostic procedures

Release

COA-0075 06 Page 1 of 1