

Product Number:	101473
Lot Number:	426818
Expiration:	24-Mar-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.37	7.35 - 7.45	PASS

Review/Approval

Robert Somogyi

QC Manager

Name

Title

Signature/Date

  27 MAR 2025

Quanterix Corporation

900 Middlesex Turnpike, Building 1, Billerica, MA 01821

techsupport@quanterix.com

©Quanterix Corp. Simoa® is a registered trademark of Quanterix Corp.

For research use only. Not for use in diagnostic procedures