

Product Number:	100453
Lot Number:	508606
Control Lot Number:	508606A
Expiration:	27-Sep-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	6.22	6.18 - 6.22	PASS

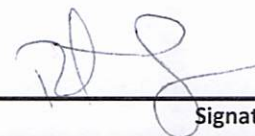
Review/Approval

Robert Somogyi

QC Manager

Name

Title



28 MAR 2021

Signature/Date

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