

<b>Product Number:</b>	100458
<b>Lot Number:</b>	502411
<b>Control Lot Number:</b>	502411A
<b>Expiration:</b>	24-Jul-2026
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Robert Somogyi

QC Manager

*Robert Somogyi* 06 MAR 2025

Name

Title

Signature/Date

Quanterix Corporation

900 Middlesex Turnpike, Building 1, Billerica, MA 01821

techsupport@quantexix.com

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

For research use only. Not for use in diagnostic procedures