

<b>Product Number:</b>	102914
<b>Lot Number:</b>	412421
<b>Expiration:</b>	3-May-2025
<b>Storage:</b>	2-8°C

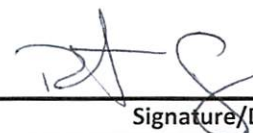
**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>	<b>PASS / FAIL</b>
pH Measurement	8.04	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi

QC Manager

 26 JUN 2024

Name

Title

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