

|                        |             |
|------------------------|-------------|
| <b>Product Number:</b> | 103847      |
| <b>Lot Number:</b>     | 313810      |
| <b>Control Number:</b> | 313810A     |
| <b>Expiration:</b>     | 18-May-2024 |
| <b>Storage:</b>        | 2 - 8°C     |

**Laboratory Analysis**

| Characteristic | Result | pH Target   |
|----------------|--------|-------------|
| pH             | Pass   | 7.35 - 7.45 |

Review/Approval

*Sonia Khillan*

QC Manager

*Sonia Khillan* 19 JUL 2023

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