

|                 |            |
|-----------------|------------|
| Product Number: | 101476     |
| Lot Number:     | 521625     |
| Expiration:     | 4-Aug-2026 |
| Storage:        | 2 - 8°C    |

Laboratory Analysis

| Characteristic | Result | pH Target   | PASS / FAIL |
|----------------|--------|-------------|-------------|
| pH             | 8.95   | 8.95 - 9.05 | PASS        |

Review/Approval

Kathleen Cornell  
Name

Quality Control  
Supervisor  
Title

Kathleen Cornell 15 Aug 2025  
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