

<b>Product Number:</b>	101956
<b>Lot Number:</b>	414207
<b>Expiration:</b>	21-Nov-2025
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.35	7.35 - 7.45	PASS

**Review/Approval**

Robert Somogyi

QC Manager

 28 MAY 2024

**Name**

**Title**

**Signature/Date**

Quanterix Corporation

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

900 Middlesex Turnpike, Building 1, Billerica, MA 01821

For research use only. Not for use in diagnostic procedures

techsupport@quanterix.com