

Product Number:	101956
Lot Number:	523313
Expiration:	21-Feb-2027
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.36	7.35 - 7.45	PASS

Review/Approval

<i>Heidi Caswell</i>	<i>Director, QC</i>	<i>Heidi Caswell</i> <i>08 Sep 2025</i>
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