

Product Number:	103208
Lot Number:	508317
Expiration:	24-Mar-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.21×10^9 Beads/mL

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title

[Signature]
10 Jun 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