

Product Number:	103208
Lot Number:	313703
Control Lot Number:	313703A
Expiration:	17-May-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.30 x 10 ⁹ Beads/mL

Review/Approval

Sonia Khillan QC Manager Sonia Khillan 17JUL2023

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