

Product Number:	103208
Lot Number:	208110
Expiration:	22-Mar-2023
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.19 x 10 <sup>9</sup> Beads/mL

Review/Approval

*Matt Smart*

Name

QC

Title

*[Signature]* 10/15/2022

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