

|                 |            |
|-----------------|------------|
| Product Number: | 103208     |
| Lot Number:     | 506320     |
| Expiration:     | 4-Mar-2026 |
| Storage:        | 2-8°C      |

**Laboratory Analysis**

| Characteristic | Means of Analysis | Result                      |
|----------------|-------------------|-----------------------------|
| Concentration  | Multisizer        | $1.32 \times 10^9$ Beads/mL |

Review/Approval

Kathleen Cornell  
Name

Quality Control  
Supervisor  
Title

 27 May 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