

Product Number:	100993
Lot Number:	512902
Expiration:	9-May-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	5.27×10^6 Beads/mL

Review/Approval

Kathleen Cornell

Name

*Quality Control
Supervisor*

Title

[Signature]
15 May 2025

Signature/Date