

Product Number:	103527
Lot Number:	410217
Expiration:	11-Apr-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.34×10^9 Beads/mL

Review/Approval

Kathleen Cornell
Name

*Quality Control
Supervisor*
Title

[Signature]
04 Dec 2025
Signature/Date