

Product Number:	103526
Lot Number:	403315
Expiration:	2-Feb-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.47×10^9 Beads/mL

Review/Approval

Kathleen Cornell

Name

Quality Control
Supervisor

Title

[Signature]
18 Aug 2025

Signature/Date