

Product Number:	103526
Lot Number:	403317
Expiration:	2-Feb-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.18×10^9 Beads/mL

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title

[Signature]
04 Dec 2025
Signature/Date