

Product Number:	103527
Lot Number:	410216
Expiration:	11-Apr-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.29×10^9 Beads/mL

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title

[Signature]
19 Nov 2025
Signature/Date