

Product Number:	103527
Lot Number:	410214
Expiration:	11-Apr-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.33×10^9 Beads/mL

Review/Approval

Kathleen
Cornell

Name

Quality Control
Supervisor

Title


10 Jun 2025

Signature/Date