

Product Number:	103208
Lot Number:	521015
Expiration:	29-Jul-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


23 Sep 2025

Signature/Date