

Product Number:	103208
Lot Number:	532308
Expiration:	19-Nov-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.20×10^9 Beads/mL

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title


04 Dec 2025
Signature/Date