

Product Number:	103208
Lot Number:	517405
Expiration:	23-Jun-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.22×10^9 Beads/mL

Review/Approval

Kathleen Cornell

Name

Quality Control
Supervisor

Title

Kathleen Cornell
30 Jul 2025

Signature/Date