

Product Number:	103208
Lot Number:	521019
Expiration:	29-Jul-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.28×10^9 Beads/mL

Review/Approval

Kathleen Cornell

Name

*Quality Control
Supervisor*

Title

[Signature]
20 Nov 2025

Signature/Date