

Product Number:	103529
Lot Number:	410213
Expiration:	11-Apr-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.13 x 10 ⁹ Beads/mL

Review/Approval

Heidi Caswell

Director, QC

H. Caswell 11 Feb 2025

Name

Title

Signature/Date