

Product Number:	103526
Lot Number:	335008
Expiration:	16-Dec-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.31×10^9 Beads/mL

Review/Approval

Robert Somogyi

QC Manager

 21 FEB 2025

Name

Title

Signature/Date