

Product Number:	103526
Lot Number:	335213
Expiration:	18-Dec-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.21 x 10 ⁹ Beads/mL

Review/Approval

ROBERT SOMOBYI
Name

QC MANAGER
Title

 24 JAN 2024
Signature/Date