

Product Number:	103526
Lot Number:	335006
Expiration:	16-Dec-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.29 x 10 ⁹ Beads/mL

Review/Approval

Robert Somogyi
Name

Manager, Quality Control
Title

 09 DEC 2024
Signature/Date