

Product Number:	103527
Lot Number:	400919
Expiration:	9-Jan-2026
Storage:	2 - 8°C


Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.16 x 10 ⁹ Beads/mL

Review/Approval

Robert Somogyi
Name

QC Manager
Title

 24SEP2024
Signature/Date