

Product Number:	103529
Lot Number:	405411
Expiration:	23-Feb-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.39 x 10 ⁹ Beads/mL

Review/Approval

Robert Somogyi

QC Manager

 27 DEC 2024

Name

Title

Signature/Date