

Product Number:	103528
Lot Number:	403314
Expiration:	2-Feb-2026
Storage:	2 - 8°C


Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.15 x 10 ⁹ Beads/mL

Review/Approval

Robert Somogyi

QC MANAGER

 01 DEC 2024

Name

Title

Signature/Date