

Product Number:	103612
Lot Number:	102535
Expiration:	5-Jun-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.39 x 10 ⁹ Beads/mL

Review/Approval

Sonia Khillan QC Manager Sonia Khillan / 01NOV2022

Name

Title

Signature/Date