

Product Number:	103526
Lot Number:	302011
Expiration:	20-Jan-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.36 x 10 ⁹ Beads/mL

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia 12/11/2023

Signature/Date