

Product Number:	103526
Lot Number:	212207
Expiration:	2-May-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.25×10^9 Beads/mL

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 17 JAN 2023
Signature/Date