

Product Number:	103526
Lot Number:	233613
Expiration:	2-Dec-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.21 x 10 ⁹ Beads/mL

Review/Approval

Sonia Khillan
Name

QC manager
Title

Soni 12/16/22 DEC 2022
Signature/Date