

Product Number:	103526
Lot Number:	015312
Expiration:	01-Jun-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.19x10 ⁹ Beads/mL

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 24Jul2020

Name

Title

Signature/Date