

Product Number:	103526
Lot Number:	026516
Expiration:	21-Sep-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.29x10 ⁹ Beads/mL

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 06Oct2021

Name

Title

Signature/Date