

Product Number:	103526
Lot Number:	124615
Expiration:	3-Sep-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.33x10 ⁹ Beads/mL

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 18Oct2021

Signature/Date