

Product Number:	103526
Lot Number:	115513
Expiration:	4-Jun-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.16x10 ⁹ Beads/mL

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 27 Jul 2021

Name

Title

Signature/Date