

Product Number:	103526
Lot Number:	123212
Expiration:	20-Aug-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.361x10 ⁹ Beads/mL

Review/Approval

Shivam Goel

Name

QC

Title

Shivam Goel 07 Dec 2021

Signature/Date