

Product Number:	103526
Lot Number:	110709
Expiration:	17-Apr-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.23x10 ⁹ Beads/mL

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 28 Apr 2021
Signature/Date