

Product Number:	103526
Lot Number:	210511
Expiration:	15-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.23×10^9 Beads/mL

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 05 May 2022
Signature/Date