

Product Number:	103526
Lot Number:	201410
Expiration:	14-Jan-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.033 x 10 ⁹ Beads/mL

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 26 Jan 2022

Signature/Date