

Product Number:	103528
Lot Number:	007218
Expiration:	12-Mar-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.26x10 ⁹ Beads/mL

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 04/20/2020

Name

Title

Signature/Date