

Product Number:	103529
Lot Number:	007219
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 09Jun2020

Name

Title

Signature/Date