

Product Number:	103529
Lot Number:	007215
Expiration:	12-Mar-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.17x10 ⁹ Beads/mL

Review/Approval

Shivani Goel

Name

QC Manager

Title

Shivani Goel 01May2020

Signature/Date