

Product Number:	103528
Lot Number:	100819
Expiration:	8-Jan-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.00x10 ⁹ Beads/mL

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 05Feb2021

Name

Title

Signature/Date