

Product Number:	103528
Lot Number:	118312
Expiration:	2-Jul-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.03x10 ⁹ Beads/mL

Review/Approval

Shwani Goel

Name

QC

Title

Shwani Goel 23 Sep 2021

Signature/Date