

Product Number:	103529
Lot Number:	201411
Expiration:	14-Jan-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.067×10^9 Beads/mL

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 27Jan2022

Signature/Date