

Product Number:	103528
Lot Number:	126312
Expiration:	20-Sep-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.34 x10 ⁹ Beads/mL

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 18Jan2022

Signature/Date