

Product Number:	101761
Lot Number:	936023
Expiration:	26-Dec-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.24x10 ⁹ Beads/mL

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 14Apr2020
Signature/Date