

Product Number:	101986
Lot Number:	931124
Expiration:	7-Nov-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.25x10 ⁹ Beads/mL

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 14Apr2020
Signature/Date