

Product Number:	101985
Lot Number:	931120
Expiration:	7-Nov-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.16x10 ⁹ Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

14 Feb 2020

Name

Title

Signature/Date