

Product Number:	101985
Lot Number:	936017
Expiration:	26-Dec-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.32x10 ⁹ Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

p.stadler 13Jan2020

Name

Title

Signature/Date