

Product Number:	101987
Lot Number:	905914
Expiration:	28-Aug-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.38x10 ⁹ Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

p.stadler 20 Feb 2020

Name

Title

Signature/Date