

Product Number:	101986
Lot Number:	005114
Expiration:	20-Feb-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.21×10^9 Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

p. stadler

28 Feb 2020

Name

Title

Signature/Date