

Product Number:	101986
Lot Number:	900713
Expiration:	7-Jul-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Bead Concentration (beads/mL)	Multisizer	1.37×10^9

Review/Approval

Patrizia stadler

QC Assoc Director

P. Stadler 27 Feb 2019

Name

Title

Signature/Date