

Product Number:	101986
Lot Number:	906017
Expiration:	1-Sep-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Bead Concentration (beads/mL)	Multisizer	1.63 x 10 ⁹

Review/Approval

Patrizia Stadler

Assoc. Director, QC

Name

Title

Signature/Date