

Product Number:	101986
Lot Number:	830514
Expiration:	1-May-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Bead Concentration (beads/mL)	Multisizer	1.28 x 10 ⁹

Review/Approval

patrizia Stadler QC Assoc. Director

Name

Title

P. Stadler 14 Feb 2019

Signature/Date