

Product Number:	101985
Lot Number:	905210
Expiration:	21-Aug-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.43×10^9 Beads/mL

Review/Approval

Patrizia Stadler

QC Assoc Director

P. Stadler 28 Feb 2019

Name

Title

Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.