

Product Number:	101761
Lot Number:	921407
Expiration:	2-Feb-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.28×10^9 Beads/mL

Review/Approval

Patrizia Stadler

QC Assoc Director

P. Stadler

25 Oct 2019

Name

Title

Signature/Date

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