

Product Number:	101987
Lot Number:	931114
Expiration:	7-May-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.18 x 10 ⁹ Beads/mL

Review/Approval Patrizia Stadler Assoc. Director, QC *P. Stadler* 15 Nov 2019
Name Title Signature/Date

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