

Product Number:	101987
Lot Number:	905915
Expiration:	28-Feb-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.37 x 10 ⁹ Beads/mL

Review/Approval Patrizia Stadler Assoc. Director, QC P. Stadler 20 Dec 2019
Name Title Signature/Date

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