

Product Number:	101986
Lot Number:	920716
Expiration:	26-Jan-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.25 x 10 ⁹ Beads/mL

Review/Approval Patrizia Stadler Assoc. Director, QC P. Stadler 05 Aug 2019
Name Title Signature/Date

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