

Product Number:	101986
Lot Number:	931116
Expiration:	7-May-2020
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.12 x 10 <sup>9</sup> 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.