

Product Number:	101987
Lot Number:	920717
Expiration:	26-Jan-2020
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.28 x 10 <sup>9</sup> 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.