

Product Number:	101986
Lot Number:	936014
Expiration:	26-Dec-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.30 x 10 ⁹ Beads/mL

Review/Approval Patrizia Stadler Assoc. Director, QC P. Stadler 07 Jan 2020
 Name Title Signature/Date

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