

Product Number:	103208
Lot Number:	915810
Expiration:	10-Dec-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.66x10 ⁹ Beads/mL

Review/Approval Patrizia Stadler Assoc. Director, QC P. Stadler 26 JUN 2019
Name Title Signature/Date

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