

Product Number:	103208
Lot Number:	926315
Expiration:	20-Mar-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.22×10^9 Beads/mL

Review/Approval Patrizia Stadler Assoc. Director, QC P. Stadler 27 Sep 2019
Name Title Signature/Date

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