

Product Number:	103208
Lot Number:	922008
Expiration:	8-Feb-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.25x10 ⁹ Beads/mL

Review/Approval Patrizia Stadler Assoc. Director, QC *P. Stadler* 09 Aug 2019
Name Title Signature/Date

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