

Product Number:	103207
Lot Number:	823621
Expiration:	30-Jan-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.44×10^9 Beads/mL

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler

27 Sep 2019

Signature/Date

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