

Product Number:	103207
Lot Number:	912104
Expiration:	30-Jan-2021
Storage Conditions:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.20×10^9 Beads/mL

Review/Approval

Patrizia Stadler

QC Assoc. Director

P. Stadler

28 May 2019

Name

Title

Signature/Date

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