

Product Number:	101985
Lot Number:	936013
Expiration:	26-Dec-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.26x10 ⁹ Beads/mL

Review/Approval

Patrizia Stadler
Name

QC Assoc Director
Title

P. Stadler

07 Jan 2020

Signature/Date

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