

Product Number:	101987
Lot Number:	936015
Expiration:	26-Dec-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.23 x 10 ⁹ Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

07 Jan 2020

Name

Title

Signature/Date

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