

Product Number:	101985
Lot Number:	005219
Expiration:	21-Feb-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.41x 10 ⁹ Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

28 Feb 2020

Name

Title

Signature/Date

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