

Product Number:	103208
Lot Number:	029314
Expiration:	19-Oct-2021
Storage:	2-8 °C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.40x 10 ⁹ Beads/mL

Review/Approval

Patrizia Stadler

QC Assoc Director

Shavani Goel 19 Nov 2020

Name

Title

Signature/Date

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