

Product Number:	103527
Lot Number:	019025
Expiration:	8-Jul-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.24 x 10 ⁹ Beads/mL

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 13 Aug 2020

Name

Title

Signature/Date

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