

Product Number:	103208
Lot Number:	026914
Expiration:	25-Sep-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.29x10 ⁹ Beads/mL

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 20 Oct 2020

Name

Title

Signature/Date

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