

Product Number:	100995
Lot Number:	018903
Expiration:	7-Jul-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	9.80x10 ⁶ Beads/mL

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shwam Goel 28 Jul 2020
Signature/Date

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