

Product Number:	103208
Lot Number:	105015
Expiration:	19-Feb-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.20×10^9 Beads/mL

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 12 Mar 2021
Signature/Date

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