

Product Number:	103611
Lot Number:	005515
Expiration:	5-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.07×10^9 Beads/mL

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 16 Apr 2020
Signature/Date

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