

Product Number:	103611
Lot Number:	005517
Expiration:	5-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	2.82x10 ⁹ Beads/mL

Review/Approval _____
 Name: Shivani Goel Title: QC Manager Signature/Date: *Shivani Goel* *sg 11 May 2020*
11 May 2020
sg 11 May 2020

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