

Product Number:	103611
Lot Number:	015311
Expiration:	5-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.30×10^9 Beads/mL

Review/Approval Shivani Goel QC Manager Shivani Goel 09 Jul 2020
 Name Title Signature/Date

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