

Product Number:	103612
Lot Number:	005518
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 02 Jun 2020.

Name

Title

Signature/Date

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