

Product Number:	103612
Lot Number:	015313
Expiration:	5-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.24×10^9 Beads/mL

Review/Approval

Shivani Goel

QC Manager

Name

Title

Shivani Goel 03 Sep 2020
Signature/Date

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