

Product Number:	103611
Lot Number:	108830
Expiration:	5-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	2.60×10^9 Beads/mL

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 15 Jun 2021
Signature/Date

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