

Product Number:	104007
Lot Number:	126705
Expiration:	24-Sep-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.53 x 10 ⁹ Beads/mL

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 15 Feb 2022
Signature/Date

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