

Product Number:	104006
Lot Number:	132310
Expiration:	19-Nov-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.35×10^9 Beads/mL

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 28 Feb 2022

Name

Title

Signature/Date

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