

Product Number:	100993
Lot Number:	205513
Expiration:	24-Feb-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	5.2 x 10 <sup>6</sup> Beads/mL

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel* 01 Mar 2022

Name

Title

Signature/Date

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