

Product Number:	103208
Lot Number:	134926
Expiration:	15-Dec-2022
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.21 x 10 <sup>9</sup> Beads/mL

Review/Approval Shivani Goel QC, Manager Shivani Goel 07Feb2022  
 Name Title Signature/Date

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