

<b>Product Number:</b>	103612
<b>Lot Number:</b>	130718
<b>Expiration:</b>	5-Dec-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.12 x10 <sup>9</sup> Beads/mL

Review/Approval

*Shivani Goel*  
Name

QC

Title

*Shivani Goel 20 Dec 2021*

Signature/Date

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