

<b>Product Number:</b>	103208
<b>Lot Number:</b>	120412
<b>Expiration:</b>	23-Jul-2022
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	1.29 x 10 <sup>9</sup> Beads/mL

Review/Approval

Shivani Goel  
Name

QC  
Title

Shivani Goel 12Oct2021  
Signature/Date

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