

<b>Product Number:</b>	103208
<b>Lot Number:</b>	021307
<b>Expiration:</b>	31-Jul-2021
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	1.04x10 <sup>9</sup> Beads/mL

**Review/Approval**

Shivani Goel

QC Manager

*Shivani Goel 01sep2020*

**Name**

**Title**

**Signature/Date**

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