

<b>Product Number:</b>	103527
<b>Lot Number:</b>	007216
<b>Expiration:</b>	12-Mar-2021
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	1.26 x 10 <sup>9</sup> Beads/mL

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel 13 May 2020*

Name

Title

Signature/Date

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