

<b>Product Number:</b>	103208
<b>Lot Number:</b>	010717
<b>Expiration:</b>	16-Oct-2020
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	$1.29 \times 10^9$ Beads/mL

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel* 09/16/2020

Name

Title

Signature/Date

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