

<b>Product Number:</b>	103611
<b>Lot Number:</b>	024812
<b>Expiration:</b>	5-Dec-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	$3.17 \times 10^9$ Beads/mL

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 10 Sep 2020  
Signature/Date

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