

<b>Product Number:</b>	101985
<b>Lot Number:</b>	931123
<b>Expiration:</b>	7-Nov-2020
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	$1.76 \times 10^9$ Beads/mL

Review/Approval

*Shivani Goel*

Name

*QC Manager*

Title

*Shivani Goel 14Apr2020*

Signature/Date