

<b>Product Number:</b>	101987
<b>Lot Number:</b>	931125
<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	1.20x10 <sup>9</sup> Beads/mL

Review/Approval

*Shwani Goel*  
Name

*QC Manager*  
Title

*Shwani Goel 14Apr2020*  
Signature/Date