

<b>Product Number:</b>	103527
<b>Lot Number:</b>	322621
<b>Expiration:</b>	14-Aug-2024
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	1.306 x 10 <sup>9</sup> Beads/mL

Review/Approval

*Pratish SURETY*  
Name

*QC SUPERVISOR*  
Title

*Pratish SURETY* 30 AUG 2023  
Signature/Date