

<b>Product Number:</b>	103528
<b>Lot Number:</b>	114118
<b>Expiration:</b>	21-May-2022
<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

Shivani Goel  
Name

QC  
Title

Shivani Goel 21 Sep 2021  
Signature/Date