

<b>Product Number:</b>	103528
<b>Lot Number:</b>	007214
<b>Expiration:</b>	12-Mar-2021
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.20x10 <sup>9</sup> Beads/mL

Review/Approval

*Shivani Goel*

*QC Manager*

*Shivani Goel 01 May 2021*

Name

Title

Signature/Date