

<b>Product Number:</b>	103529
<b>Lot Number:</b>	112011
<b>Expiration:</b>	30-Apr-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.38x10 <sup>9</sup> Beads/mL

Review/Approval

Shivani Goel  
Name

QC

Title

Shivani Goel 07 Jul 2021  
Signature/Date