

<b>Product Number:</b>	103526
<b>Lot Number:</b>	104915
<b>Expiration:</b>	18-Feb-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.16x10 <sup>9</sup> Beads/mL

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 26Mar2021  
Signature/Date