

<b>Product Number:</b>	100998
<b>Lot Number:</b>	020208
<b>Expiration:</b>	20-Jul-2021
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	5.12x10 <sup>6</sup> Beads/mL

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel 27 Jul 2020*

Name

Title

Signature/Date