

<b>Product Number:</b>	100993
<b>Lot Number:</b>	319409
<b>Control Number:</b>	319409A
<b>Expiration:</b>	13-Jul-2024
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	4.98 x 10 <sup>6</sup> Beads/mL

Review/Approval

Sonia Khillan  
Name

QC Manager  
Title

Soni K/F/L 20 JUL 2023  
Signature/Date