

<b>Product Number:</b>	100993
<b>Lot Number:</b>	415008
<b>Expiration:</b>	29-May-2025
<b>Storage:</b>	2 - 8°C


**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	4.99 x 10 <sup>6</sup> Beads/mL

Review/Approval

Robert Somogyi  
Name

QC Manager  
Title

 13 JUN 2024  
Signature/Date