

<b>Product Number:</b>	103528
<b>Lot Number:</b>	231419
<b>Expiration:</b>	10-Nov-2023
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	1.40 x 10 <sup>9</sup> Beads/mL

Review/Approval

Sonia Khillan  
Name

QC Manager  
Title

Sonia R. K. / 11 JAN 2023  
Signature/Date