

<b>Product Number:</b>	103527
<b>Lot Number:</b>	210512
<b>Expiration:</b>	15-Apr-2023
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	1.48 x 10 <sup>9</sup> Beads/mL

Review/Approval

Matt Smart  
Name

QC  
Title

MSA 19 May 2022  
Signature/Date