

<b>Product Number:</b>	103529
<b>Lot Number:</b>	112010
<b>Expiration:</b>	30-Apr-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	$1.34 \times 10^9$ Beads/mL

Review/Approval

Shivani Goel  
Name

QC  
Title

Shivani Goel 27 May 2021  
Signature/Date