

<b>Product Number:</b>	103528
<b>Lot Number:</b>	209410
<b>Expiration:</b>	4-Apr-2023
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	1.32 x10 <sup>9</sup> Beads/mL

[Empty rectangular box for additional notes or comments]

Review/Approval

Richa Shrestha  
Name

QC Supervisor  
Title

[Signature] 26 Apr 2022  
Signature/Date