

<b>Product Number:</b>	103526
<b>Lot Number:</b>	120033
<b>Expiration:</b>	19-Jul-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	1.22x10 <sup>9</sup> Beads/mL

Review/Approval

*Jungil Park*

Name

*QC supervisor*

Title

*[Signature]* 01/29/2021

Signature/Date