

<b>Product Number:</b>	103208
<b>Lot Number:</b>	221512
<b>Expiration:</b>	3-Aug-2023
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.21 x 10 <sup>9</sup> Beads/mL

Review/Approval

*Sonia Khillan*  
Name

*QC Manager*  
Title

*Sonia Khillan / 28 DEC 2022*  
Signature/Date