

<b>Product Number:</b>	103528
<b>Lot Number:</b>	304809
<b>Expiration:</b>	17-Feb-2024
<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

*Pratish Seneviratne*  
Name

*QC*  
*SUPERVISOR*  
Title

*Pratish Seneviratne*  
Signature/Date  
*12 OCT 2022*