

Product Number:	103529
Lot Number:	324410
Expiration:	1-Sep-2024
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.30 x 10 <sup>9</sup> Beads/mL

Review/Approval

*Sonia Khillan*  
Name

*QC Manager*  
Title

*Sonia Khillan 13 OCT 2023*  
Signature/Date