

Product Number:	103529
Lot Number:	119716
Expiration:	16-Jul-2022
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.41 X 10 <sup>9</sup> Beads/mL

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel* 04 Apr 2022

Name

Title

Signature/Date