

Product Number:	103208
Lot Number:	317112
Control Number:	317112A
Expiration:	20-Jun-2024
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.18 x 10 <sup>9</sup> Beads/mL

Review/Approval

Sonia Khillan                      QC Manager                      Sonia Khillan 23 Aug 2023  
Name    Title    Signature/Date