

Product Number:	103208
Lot Number:	406010
Expiration:	28-Feb-2025
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.23 x 10 <sup>9</sup> Beads/mL

Review/Approval

Robert Somogyi  
Name

QC Manager  
Title

 09 JUL 2024  
Signature/Date