

Product Number:	103208
Lot Number:	505812
Expiration:	27-Feb-2026
Storage:	2-8°C

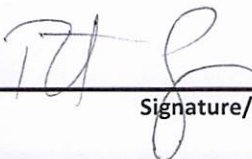
**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.19 x 10 <sup>9</sup> Beads/mL

Review/Approval

Robert Somogyi  
Name

QC Manager  
Title

 11 MAR 2025  
Signature/Date