

Product Number:	103208
Lot Number:	506315
Expiration:	4-Mar-2026
Storage:	2-8°C


**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	$1.32 \times 10^9$ Beads/mL

Review/Approval

Robert Somogyi  
Name

QC Manager  
Title

  
Signature/Date

24 MAR 2025