

Product Number:	103208
Lot Number:	308308
Expiration:	24-Mar-2024
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	$1.35 \times 10^9$ Beads/mL

Review/Approval

Sonia Khillan  
Name

QC Manager  
Title

Sonia Khillan 22 JUN 2023  
Signature/Date