

<b>Product Number:</b>	103527
<b>Lot Number:</b>	302705
<b>Expiration:</b>	27-Jan-2024
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	$1.56 \times 10^9$ Beads/mL

Review/Approval

Sonia Khilkin  
Name

QC Manager  
Title

[Signature] 11APR2023  
Signature/Date