

<b>Product Number:</b>	101987
<b>Lot Number:</b>	936019
<b>Expiration:</b>	26-Dec-2020
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	$1.24 \times 10^9$ Beads/mL

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*P. Stadler* 13 Jan 2020  
Signature/Date