

<b>Product Number:</b>	101985
<b>Lot Number:</b>	905213
<b>Expiration:</b>	21-Feb-2020
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	1.45x10 <sup>9</sup> Beads/mL

**Review/Approval**

Patrizia Stadler  
**Name**

Assoc. Director, QC  
**Title**

*P Stadler 20Jan2020*  
**Signature/Date**