

<b>Product Number:</b>	101761
<b>Lot Number:</b>	936021
<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	1.22x10 <sup>9</sup> Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

*p. Stadler* 14 Feb 2020

Name

Title

Signature/Date