

<b>Product Number:</b>	101986
<b>Lot Number:</b>	936018
<b>Expiration:</b>	26-Dec-2020
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.24x10 <sup>9</sup> Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

*P. Stadler 13 Jan 2020*

Name

Title

Signature/Date