

<b>Product Number:</b>	101761
<b>Lot Number:</b>	921409
<b>Expiration:</b>	2-Aug-2020
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	$1.28 \times 10^9$ Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

*P. Stadler* 28 Feb 2020

Name

Title

Signature/Date