

<b>Product Number:</b>	101761
<b>Lot Number:</b>	906515
<b>Expiration:</b>	6-Sep-2019
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	1.34x10 <sup>9</sup> Beads/mL

Review/Approval

Patrizia Stadler

QC Assoc Director

*P. Stadler* 18 Jul 2019

Name

Title

Signature/Date

**Statement:** The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.