

<b>Product Number:</b>	101761
<b>Lot Number:</b>	920605
<b>Expiration:</b>	25-Jan-2020
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	1.36x10 <sup>9</sup> Beads/mL

Review/Approval

Patrizia Stadler

QC Assoc Director

*P. Stadler*

09 Oct 2019

Name

Title

Signature/Date

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