

<b>Product Number:</b>	101987
<b>Lot Number:</b>	901811
<b>Expiration:</b>	18-Jul-2019
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Bead Concentration (beads/mL)	Multisizer	1.18 x 10 <sup>9</sup>

Review/Approval

Patrizia Stadler QC Assoc Director

*P. Stadler* 21 Feb 2019

Name

Title

Signature/Date

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