

<b>Product Number:</b>	101987
<b>Lot Number:</b>	905911
<b>Expiration:</b>	28-Aug-2019
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Bead Concentration (beads/mL)	Multisizer	1.49 x 10 <sup>9</sup>

**Review/Approval**

Patrizia Stadler

QC Assoc Director

*P. Stadler* 07 Mar 2019

**Name**

**Title**

**Signature/Date**

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