

<b>Product Number:</b>	101987
<b>Lot Number:</b>	817008
<b>Expiration:</b>	19-Dec-2018
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Bead Concentration (beads/mL)	Multisizer	$1.29 \times 10^9$

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

 05 Dec 2018

Signature/Date

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