


<b>Product Number:</b>	101761
<b>Lot Number:</b>	819910
<b>Expiration:</b>	18-Jan-2019
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	$1.20 \times 10^9$ Beads/mL

<b>Review/Approval</b>	<u>Sivasankari Murugan</u>	<u>QC Supervisor</u>	<u> 05 OCT 2018</u>
	<b>Name</b>	<b>Title</b>	<b>Signature/Date</b>

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