

<b>Product Number:</b>	103208
<b>Lot Number:</b>	927412
<b>Expiration:</b>	1-Apr-2020
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	$1.31 \times 10^9$ Beads/mL

Review/Approval      Patrizia Stadler      Assoc. Director, QC      P. Stadler      25 Oct 2019  
Name      Title      Signature/Date

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