

<b>Product Number:</b>	103208
<b>Lot Number:</b>	936104
<b>Expiration:</b>	27-Jun-2020
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	1.35x10 <sup>9</sup> Beads/mL

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*P. Stadler*

31 Jan 2020

Signature/Date

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