

<b>Product Number:</b>	103208
<b>Lot Number:</b>	922008
<b>Expiration:</b>	8-Feb-2020
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.25x10 <sup>9</sup> Beads/mL

Review/Approval Patrizia Stadler Assoc. Director, QC P. Stadler 09 Aug 2019  
Name Title Signature/Date

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