

<b>Product Number:</b>	101985
<b>Lot Number:</b>	905212
<b>Expiration:</b>	21-Feb-2020
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.39x10 <sup>9</sup> Beads/mL

<b>Review/Approval</b>	Patrizia Stadler	QC Assoc Director	<i>P. Stadler</i>	20 Dec 2019
	Name	Title	Signature/Date	

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