

<b>Product Number:</b>	100993
<b>Lot Number:</b>	010007
<b>Expiration:</b>	9-Apr-2021
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	$4.78 \times 10^6$ Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

*P. Stadler*

14 Apr 2020

Name

Title

Signature/Date

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