

Product Number:	103206
Lot Number:	823620
Expiration:	30-Jan-2021
Storage Conditions:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	$3.41 \times 10^9$ Beads/mL

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*P. Stadler* 29 Nov 2019  
Signature/Date

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