

Product Number:	103206
Lot Number:	823622
Expiration:	30-Jan-2021
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	$3.62 \times 10^9$ Beads/mL

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*P. Stadler*

18 Oct 2019

Signature/Date

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