

Product Number:	103317
Lot Number:	834713
Expiration:	27-Nov-2019
Storage Conditions:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	2.67×10^9 Beads/mL

Review/Approval

Patrizia Stadler

Associate Director, QC

Name

Title

P. Stadler 1/9/2018
Signature/Date

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