

Product Number:	103208
Lot Number:	829017
Expiration:	17-Apr-2019
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.37×10^9 Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler 27 Feb 2019

Name

Title

Signature/Date

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