

Product Number:	103208
Lot Number:	907111
Expiration:	12-Sep-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.45×10^9 Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

Name

Title

P. Stadler

25 JUN 2019

Signature/Date

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