

Product Number:	103208
Lot Number:	928412
Expiration:	11-Apr-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.25×10^9 Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

25NOV2019

Name

Title

Signature/Date

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