

Product Number:	103206
Lot Number:	001510
Expiration:	30-Jan-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.24×10^9 Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

22 Jan 2020

Name

Title

Signature/Date

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