

Product Number:	101761
Lot Number:	906509
Expiration:	6-Sep-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.30×10^9 Beads/mL

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler 12/10/2019

Name

Title

Signature/Date

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