

Product Number:	101985
Lot Number:	831905
Expiration:	15-May-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.35×10^9 Beads/mL

Review/Approval Patrizia Stadler QC Assoc Director p. Stadler 21 Feb 2019
Name Title Signature/Date

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