

Product Number:	101985
Lot Number:	905211
Expiration:	21-Aug-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.41x10 ⁹ Beads/mL

Review/Approval	Patrizia Stadler Name	Assoc. Director, QC Title	<i>P. Stadler</i> Signature	03 JUL 2019 Date
-----------------	--------------------------	------------------------------	--------------------------------	---------------------

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