

Product Number:	103208
Lot Number:	923206
Expiration:	20-Feb-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.31x10 ⁹ Beads/mL

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	05 Sep 2019
	Name	Title	Signature/Date	

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