

Product Number:	103208
Lot Number:	903908
Expiration:	8-Aug-2019
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.42x10 ⁹ Beads/mL

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i> 27 Feb 2019
	<small>Name</small>	<small>Title</small>	<small>Signature/Date</small>

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