

Product Number:	101761
Lot Number:	920606
Expiration:	25-Jul-2020
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.36x10 ⁹ Beads/mL

Review/Approval	Patrizia Stadler	QC Assoc Director	<i>P. Stadler</i>	20 Dec 2019
	Name	Title	Signature/Date	

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