

Product Number:	103208
Lot Number:	007307
Expiration:	13-Sep-2020
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.32x10 ⁹ Beads/mL

Review/Approval

	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	25 Mar 2020
	Name	Title	Signature	Date

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