

Product Number:	103527
Lot Number:	019027
Expiration:	8-Jul-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.35 x 10 ⁹ Beads/mL

Review/Approval	Shivani Goel	QC Manager	<i>Shivani Goel</i> 19 Nov 2020
	Name	Title	Signature/Date

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