

Product Number:	103612
Lot Number:	024813
Expiration:	5-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.28×10^9 Beads/mL

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 200ct2020
	Name	Title	Signature/Date

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