

Product Number:	103611
Lot Number:	024814
Expiration:	5-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	2.82x10 ⁹ Beads/mL

Review/Approval	Shivani Goel Name	QC Manager Title	<i>Shivani Goel 20 Oct 2020</i> Signature/Date
-----------------	----------------------	---------------------	---

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