

Product Number:	103208
Lot Number:	019111
Expiration:	9-Jan-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.25x10 ⁹ Beads/mL

Review/Approval

Shivani Goel	QC Manager	<i>Shivani Goel</i> 20 Jul 2020
Name	Title	Signature/Date

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