

Product Number:	104006
Lot Number:	206617
Expiration:	7-Mar-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.18 x 10 ⁹ Beads/mL

Note: Beads provided are conjugated with 488 nm fluorescent dye.
For more information, please visit
<https://go.quanterix.com/HomebrewBeadUpdate>



Review/Approval _____
Name: Shivani Goel Title: QC Manager Signature/Date: *Shivani Goel 23 Mar 2022*

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