

Product Number:	103612
Lot Number:	130715
Expiration:	5-Dec-2022
Storage:	2 - 8 °C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	2.79 x10 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

Shivani Goel
Name

QC

Title

Shivani Goel 29 NOV 2021
Signature/Date

Signature/Date

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