

Product Number:	104006
Lot Number:	131517
Expiration:	11-Nov-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.33 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 28 Dec 2021
Signature/Date

Signature/Date

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