

Product Number:	104006
Lot Number:	131609
Expiration:	12-Nov-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.31 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 Gael
Name

QC

Title

Shivani Gael 28 Dec 2021

Signature/Date

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