

Product Number:	104006
Lot Number:	312409
Control Number:	312409A
Expiration:	4-May-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.30×10^9 Beads/mL

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



Review/Approval

Pratish SNETI

Name

QC SUPERVISOR

Title

Pratish SNETI

Signature/Date

09 AUG 2023