

Product Number:	104006
Lot Number:	308219
Expiration:	23-Mar-2025
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

Brian Ciccarelli

Name

VP, QA + RA

Title

B-C

Signature/Date

19 APR 2023