

Product Number:	104007
Lot Number:	312408
Control Number:	312408A
Expiration:	4-May-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.27 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 Sonia Khillan QC Manager Sonia Khillan 27 JUL 2023

Name Title Signature/Date