

Product Number:	104006
Lot Number:	304013
Expiration:	9-Feb-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.33 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

Name

QC Manager

Title

Soni 12/11. 17.MAR2023

Signature/Date