

Product Number:	104006
Lot Number:	231817
Expiration:	14-Nov-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.25 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 Khillar

Name

QC Manager

Title

Sonia Khillar 13 DEC 2022

Signature/Date