

Product Number:	104007
Lot Number:	303713
Expiration:	6-Feb-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.23 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

Sania Khillan
Name

QC Manager
Title

Sania Khillan 24 FEB 2023
Signature/Date