

Product Number:	104007
Lot Number:	230417
Expiration:	31-Oct-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 Khillan QC Manager Sonia Khillan 01DEC2022

Name Title Signature/Date