

Product Number:	104007
Lot Number:	223713
Expiration:	25-Aug-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.468 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 Matt Kodouris QC Supervisor [Signature] 09Feb2023
Name
Title
Signature/Date