

Product Number:	104006
Lot Number:	301314
Expiration:	13-Jan-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.18 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 Koulavris

QC Supervisor

[Signature] 07Feb2023

Name

Title

Signature/Date