

Product Number:	104007
Lot Number:	208008
Expiration:	21-Mar-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.25 x10 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

Shivani Goel

QC Manager

Shivani Goel 20 Apr 2022

Name

Title

Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.