

Product Number:	104006
Lot Number:	206911
Expiration:	10-Mar-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.26 x 10 ⁹ Beads/mL

Note: Beads provided are conjugated with 488 nm fluorescent dye.
For more information, please visit
<https://go.quoterix.com/HomebrewBeadUpdate>



Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 04Apr2022
Signature/Date

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