

Product Number:	104006
Lot Number:	205909
Expiration:	28-Feb-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.17 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

Matthew Smart
~~Shivani Goel~~

QC Director
~~QC Manager~~

[Signature] 16Mar2022

Name *16Mar2022*

Title *16Mar2022*

Signature/Date

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