

| | |
|------------------------|------------|
| Product Number: | 103612 |
| Lot Number: | 115515 |
| Expiration: | 5-Dec-2022 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|-----------------------|--------------------------|--------------------------------|
| Concentration | Multisizer | 2.67 x10 ⁹ Beads/mL |

Review/Approval

Wenkai Tsai

Name

QC

Title



18 Aug 2021

Signature/Date

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