

|                        |            |
|------------------------|------------|
| <b>Product Number:</b> | 103611     |
| <b>Lot Number:</b>     | 130714     |
| <b>Expiration:</b>     | 5-Dec-2022 |
| <b>Storage:</b>        | 2 - 8°C    |

**Laboratory Analysis**

| Characteristic | Means of Analysis | Result                        |
|----------------|-------------------|-------------------------------|
| Concentration  | Multisizer        | 2.81x10 <sup>9</sup> Beads/mL |

Review/Approval Shivani Goel QC Shivani Goel 16 Nov 2021  
Name Title Signature/Date

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