

|                        |             |
|------------------------|-------------|
| <b>Product Number:</b> | 104007      |
| <b>Lot Number:</b>     | 233208      |
| <b>Expiration:</b>     | 28-Nov-2024 |
| <b>Storage:</b>        | 2 - 8°C     |

**Laboratory Analysis**

| Characteristic | Means of Analysis | Result                          |
|----------------|-------------------|---------------------------------|
| Concentration  | Multisizer        | 1.23 x 10 <sup>9</sup> Beads/mL |

Note: Beads provided are conjugated with 488 nm fluorescent dye.  
For more information, please visit

<https://go.quanterix.com/HomebrewBeadUpdate>



Review/Approval

*Sonia Khillan*

Name

*QC Manager*

Title

*Sonia 12/16 / 03 JAN 2023*

Signature/Date