

|                        |             |
|------------------------|-------------|
| <b>Product Number:</b> | 104007      |
| <b>Lot Number:</b>     | 307212      |
| <b>Expiration:</b>     | 13-Mar-2025 |
| <b>Storage:</b>        | 2 - 8°C     |

**Laboratory Analysis**

| <b>Characteristic</b> | <b>Means of Analysis</b> | <b>Result</b>                   |
|-----------------------|--------------------------|---------------------------------|
| Concentration         | Multisizer               | 1.32 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 QC Manager Sonia Khillan 07 APR 2023  
Name Title Signature/Date