

<b>Product Number:</b>	104006
<b>Lot Number:</b>	226507
<b>Expiration:</b>	22-Sep-2024
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
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

*Matt Koubouris*

Name

*QC Supervisor*

Title

*[Signature]* 10092622

Signature/Date