

<b>Product Number:</b>	104006
<b>Lot Number:</b>	311005
<b>Control Number:</b>	311005A
<b>Expiration:</b>	20-Apr-2025
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.27 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

*Pratish SINGH*

Name

*QC SUPERVISOR*

Title

*Pratish Singh*

Signature/Date

15 MAY 2023