

<b>Product Number:</b>	104007
<b>Lot Number:</b>	220210
<b>Expiration:</b>	21-Jul-2024
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Means of Analysis</b>	<b>Result</b>
Concentration	Multisizer	1.34 x 10 <sup>9</sup> Beads/mL

Note: Beads provided are conjugated with 488 nm fluorescent dye.  
For more information, please visit  
<https://go.quoterix.com/HomebrewBeadUpdate>



Review/Approval

*Matt Smart*  
Name

QC  
Title

*MJA 15 Aug 2022*  
Signature/Date