

<b>Product Number:</b>	104006
<b>Lot Number:</b>	228609
<b>Expiration:</b>	13-Oct-2024
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.15 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 / 01/06/2022  
 Name Title Signature/Date