

Product Number:	104006
Lot Number:	220211
Expiration:	21-Jul-2024
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
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.quanterix.com/HomebrewBeadUpdate>



Review/Approval

*Matt Smart*  
Name

QC

Title

*[Signature]* 15 Aug 2022  
Signature/Date