

Product Number:	104006
Lot Number:	231420
Expiration:	10-Nov-2024
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.22 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

*Sonia Khilkin*  
Name

*QC Manager*  
Title

*Sonia Khilkin / 11 JAN 2023*  
Signature/Date