

Product Number:	104006
Lot Number:	320907
Expiration:	28-Jul-2025
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	$1.15 \times 10^9$ 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  
Name

QC Manager  
Title

Sonia Khillan 05 SEP 2023  
Signature/Date