

Product Number:	104006
Lot Number:	330803
Expiration:	4-Nov-2025
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

**RESULT**

Multisizer Concentration	1.17 x 10 <sup>9</sup> Beads/mL
--------------------------	---------------------------------

Note: Beads provided are conjugated with 488 nm fluorescent dye.

Review/Approval

*Sonia Khillan*

Name

*Senior QC Analyst*

Title

*Sonia Khillan 01DEC2023*

Signature/Date