

Product Number:	104006
Lot Number:	418712
Expiration:	5-Jul-2026
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

**RESULT**

Multisizer Concentration	1.21 x 10 <sup>9</sup> Beads/mL
--------------------------	---------------------------------

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

Review/Approval

Robert Somogyi

Manager, Quality Control

 04DEC2024

Name

Title

Signature/Date