

Product Number:	104007
Lot Number:	534501
Expiration:	1-Dec-2027
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	$1.19 \times 10^9$ Beads/mL
--------------------------	-----------------------------

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

Review/Approval

Kathleen Corneli

Name

Quality Control  
Supervisor

Title

*[Signature]*  
15 Dec 2025

Signature/Date