

Product Number:	104006
Lot Number:	526907
Expiration:	26-Sep-2027
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	1.24×10^9 Beads/mL
--------------------------	-----------------------------

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

Review/Approval

Kathleen Cornell

Name

Quality Control
Supervisor

Title

[Signature]
21 Oct 2025

Signature/Date