

Product Number:	104006
Lot Number:	526906
Expiration:	26-Sep-2027
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	1.16×10^9 Beads/mL
--------------------------	-----------------------------

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

Review/Approval

Grey Simone

Name

Quality events and Training Specialist

Title

[Signature] 09 Oct 2025

Signature/Date