

Product Number:	104006
Lot Number:	531805
Expiration:	14-Nov-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


04 Dec 2025
Signature/Date