

Product Number:	104007
Lot Number:	519910
Expiration:	18-Jul-2027
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	1.26 x 10 ⁹ Beads/mL
--------------------------	---------------------------------

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

Review/Approval

Kathleen Cornell

Name

Quality Control
Supervisor

Title

Kathleen Cornell
15 Aug 2025

Signature/Date