

Product Number:	104006
Lot Number:	512108
Expiration:	1-May-2027
Storage:	2-8°C

Laboratory Analysis		
Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS


RESULT	
Multisizer Concentration	1.35 x 10 ⁹ Beads/mL

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

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title


24Jul2025
Signature/Date