

Product Number:	104007
Lot Number:	505210
Expiration:	21-Feb-2027
Storage:	2-8°C

Laboratory Analysis


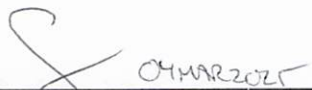
Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	1.33 x 10 ⁹ Beads/mL
--------------------------	---------------------------------

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

Review/Approval

Robert Somogyi	QC Manager		
Name	Title	Signature	Date