

Product Number:	104007
Lot Number:	503707
Control Lot Number:	503707A
Expiration:	6-Feb-2027
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT


Multisizer Concentration	1.29 x 10 ⁹ Beads/mL
--------------------------	---------------------------------

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

Review/Approval

ROBERT SOMOGYI
Name

MANAGER QUALITY CONTROL
Title

 14 FEB 2025
Signature/Date