

Product Number:	104006
Lot Number:	503706
Control Lot Number:	503706A
Expiration:	6-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

Manager, Quality Control

Name

Title

 14 FEB 2025
Signature/Date