

Product Number:	104006
Lot Number:	331510
Expiration:	11-Nov-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT


Multisizer Concentration	1.34 x 10 ⁹ Beads/mL
--------------------------	---------------------------------

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

Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

 25 JAN 2024
Signature/Date