

Product Number:	104006
Lot Number:	407413
Expiration:	14-Mar-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	1.21 x 10 ⁹ Beads/mL
--------------------------	---------------------------------

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

Review/Approval

ROBERT SOMOSYI

Name

QC MANAGER

Title

[Handwritten Signature] 16 APR 2024

Signature/Date