

Product Number:	104007
Lot Number:	334113
Expiration:	7-Dec-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	1.18 x 10 ⁹ Beads/mL
--------------------------	---------------------------------

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

Review/Approval

ROBERT SOMOSZI

Name

QC MANAGER

Title

 14 FEB 2024

Signature/Date