

Product Number:	104006
Lot Number:	333608
Expiration:	2-Dec-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

Sonia Khillan

Senior QC Analyst

Sonia Khillan 21DEC2023

Name

Title

Signature/Date