

Product Number:	104006
Lot Number:	424909
Expiration:	5-Sep-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	1.29 x 10 ⁹ Beads/mL
---------------------------------	---------------------------------

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

Review/Approval

Robert Somogyi

QC Manager

 09 Oct 2024

Name

Title

Signature/Date