

Product Number:	104006
Lot Number:	423610
Expiration:	23-Aug-2026
Storage:	2-8°C

Laboratory Analysis

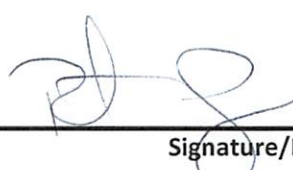
Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

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	
Name	Title	Signature/Date

27 DEC 2024