

Product Number:	104007
Lot Number:	502412
Control Lot Number:	502412A
Expiration:	24-Jan-2027
Storage:	2-8°C

Laboratory Analysis		
Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT	
Multisizer Concentration	1.22 x 10 ⁹ Beads/mL

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

Review/Approval

Robert Somogyi

QC Manager

 25 Mar 2025

Name

Title

Signature/Date