

Product Number:	104007
Lot Number:	403913
Expiration:	8-Feb-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	1.14 x 10 <sup>9</sup> Beads/mL
--------------------------	---------------------------------

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

Review/Approval

Cody Carlson  
Name

QC Manager  
Title

 07MAR2024  
Signature/Date