

Product Number:	104007
Lot Number:	513911
Expiration:	19-May-2027
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	1.20×10^9 Beads/mL
--------------------------	-----------------------------

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

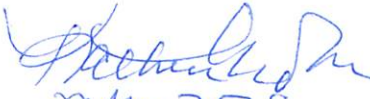
Review/Approval

Kathleen
Cornell

Name

Quality Control
Supervisor

Title


21 May 2025

Signature/Date