

Product Number:	104007
Lot Number:	507619
Expiration:	17-Mar-2027
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	1.24×10^9 Beads/mL
--------------------------	-----------------------------

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

Review/Approval

Robert Somogyi

QC Manager

Name

Title



Signature/Date