

Product Number:	104006
Lot Number:	412720
Expiration:	6-May-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

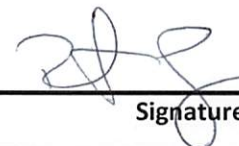
Multisizer Concentration	1.21 x 10 <sup>9</sup> Beads/mL
--------------------------	---------------------------------

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

Review/Approval

Robert Somogyi

QC Manager



28 MAY 2024

Name

Title

Signature/Date