

Product Number:	104006
Lot Number:	415218
Expiration:	31-May-2026
Storage:	2-8°C

Laboratory Analysis

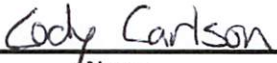
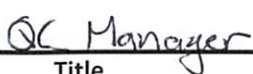

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	1.18 x 10 <sup>9</sup> Beads/mL
--------------------------	---------------------------------

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

Review/Approval

Name Title Signature/Date