

Product Number:	104006
Lot Number:	507620
Expiration:	17-Mar-2027
Storage:	2-8°C

Laboratory Analysis		
Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

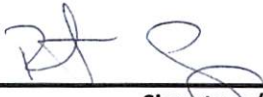
RESULT	
Multisizer Concentration	1.27 x 10 <sup>9</sup> Beads/mL

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

Review/Approval

Robert Somogyi

QC Manager

 23 APR 2025

Name

Title

Signature/Date