

Product Number:	104006
Lot Number:	522307
Expiration:	11-Aug-2027
Storage:	2-8°C

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

RESULT	
Multisizer Concentration	1.34 x 10 <sup>9</sup> Beads/mL

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

Review/Approval

Heidi Caswell

Director, QC

Heidi Caswell 08 Sep 2025

Name

Title

Signature/Date