

Product Number:	104006
Lot Number:	503708
Control Lot Number:	503708A
Expiration:	6-Feb-2027
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	1.22 x 10 <sup>9</sup> Beads/mL
--------------------------	---------------------------------

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


Review/Approval

Robert Somogyi

QC Manager

Name

Title



21 FEB 2025

Signature/Date