

Product Number:	104007
Lot Number:	329215
Expiration:	19-Oct-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	1.23 x 10 <sup>9</sup> Beads/mL
--------------------------	---------------------------------

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

Review/Approval

<u>Sonia Khillan</u>	<u>Senior QC Analyst</u>	<u>Sonia Pellera 16 NOV 2023</u>
Name	Title	Signature/Date