

Product Number:	104007
Lot Number:	223713
Expiration:	25-Aug-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

RESULT

Multisizer Concentration	1.28 x 10 ⁹ Beads/mL
--------------------------	---------------------------------

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

Review/Approval

<i>Sonia Khillan</i>	<i>Senior QC Analyst</i>	<i>Sonia Khillan 07 DEC 2023</i>
Name	Title	Signature/Date