

<b>Product Number:</b>	104006
<b>Lot Number:</b>	411620
<b>Expiration:</b>	25-Apr-2026
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	PASS / FAIL
Concentration	Multisizer	PASS

**RESULT**

<b>Multisizer Concentration</b>	1.19 x 10 <sup>9</sup> Beads/mL
---------------------------------	---------------------------------

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

Review/Approval

Robert Somogyi

QC Manager



26 JUN 2024

Name

Title

Signature/Date