

|                 |            |
|-----------------|------------|
| Product Number: | 104006     |
| Lot Number:     | 406411     |
| Expiration:     | 4-Mar-2026 |
| Storage:        | 2-8°C      |

Laboratory Analysis

| Characteristic | Means of Analysis | PASS / FAIL |
|----------------|-------------------|-------------|
| Concentration  | Multisizer        | PASS        |

RESULT

|                          |                                 |
|--------------------------|---------------------------------|
| Multisizer Concentration | 1.16 x 10 <sup>9</sup> Beads/mL |
|--------------------------|---------------------------------|

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

Review/Approval

ROBERT SOMOGYI  
Name

QC MANAGER  
Title

 15 MAR 2024  
Signature/Date