

| | |
|-----------------|-------------|
| Product Number: | 103187 |
| Lot Number: | 501524 |
| Expiration: | 25-Apr-2019 |

| Component | Item Number | Lot Number |
|----------------|-------------|------------------|
| Rainbow Beads | 101069 | MF-030518-04 |
| Low AEB Beads | 103188 | 823202 |
| High AEB Beads | 101070 | 823203 |
| RGP Reagent | 103159 | N/A ¹ |

¹ RGP is not Kit Lot Specific

Rainbow Beads

| | |
|-----------------------|------------------------|
| Bead Count (beads/mL) | 2.39 x 10 ⁶ |
| Bead % Monomeric | 91% |

Low AEB Beads

| | |
|-----------------------|------------------------|
| Bead Count (beads/mL) | 2.40 x 10 ⁶ |
| Bead % Monomeric | 90% |

High AEB Beads

| | |
|-----------------------|------------------------|
| Bead Count (beads/mL) | 2.39 x 10 ⁶ |
| Bead % Monomeric | 92% |

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title


Signature/Date