

Product Number: 101537
Lot Number: 502529
Expiration: 3-Aug-2021

| Component | Part Number | Lot Number |
|------------------|-------------|------------------|
| Bead Reagent | 101535 | 024601 |
| Detector Reagent | 101533 | 024801 |
| SBG Reagent | 101534 | 024706 |
| Control 1 | 101543 | 025205 |
| Control 2 | 101544 | 025205 |
| RGP Reagent | 103159 | N/A ¹ |

¹ RGP is not Kit Lot Specific

| | |
|-----------------------|----------------------|
| Bead Count (beads/mL) | 1.59x10 ⁷ |
| Bead % Monomeric | 93% |
| Isingle | 131 |

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 17 Sep 2020

Name

Title

Signature/Date