

| | |
|------------------------|-------------|
| Product Number: | 101537 |
| Lot Number: | 501785 |
| Expiration: | 16-Apr-2020 |

| Component | Item Number | Lot Number |
|------------------|---------------|------------------|
| Bead Reagent | 101535 | 910906 |
| Detector Reagent | 101533 | 910806 |
| SBG Reagent | 101534 | 910807 |
| Controls | 101543-101544 | 911601 |
| RGP Reagent | 103159 | N/A ¹ |

¹ RGP is not Kit Lot Specific

| | |
|-----------------------|------------------------|
| Bead Count (beads/mL) | 1.58 x 10 ⁷ |
| Bead % Monomeric | 93% |
| Isingle | 166 |

Review/Approval

Patrizia stadler QC Assoc. Director

Name

Title

p. stadler 07 MAY 2019

Signature/Date