

| | |
|-----------------|-------------|
| Product Number: | 101351 |
| Lot Number: | 502178 |
| Expiration: | 25-Aug-2020 |

| Component | Part Number | Lot Number | Concentration |
|-------------------------------|---------------|------------------|-------------------------------|
| Beads | 103206 | 823622 | 3.62X10 ⁹ beads/mL |
| Detector/Sample Diluent | 101353 | 928211 | |
| Bead Diluent | 100458 | 928215 | |
| SBG Dilution Buffer | 100376 | 930220 | |
| SBG Concentrate | 103397 | 930316 | 50nM |
| Bead Conjugation Buffer | 100453 | 932512 | |
| Bead Wash Buffer | 101352 | 928911 | |
| Bead Blocking Buffer | 100457 | 926611 | |
| Biotinylation Reaction Buffer | 100460 | 924107 | |
| EDC, 10mg | 100022 | RM2019120000025 | |
| EZ-Link NHS-PEG4 Biotin | 100348 | RM2019120000049 | |
| RGP | 103159 | N/A ¹ | |
| PSA Capture Antibody | 100463 | 930904 | 5.0 mg/mL |
| PSA Detector Antibody | 100464 | 931607 | 1.0 mg/mL |
| Calibrator A-H | 100758-100765 | 921201 | |
| Controls 1 and 2 | 100768-100769 | 921202 | |

¹ RGP is not Kit Lot specific.

| | Mean Range (pg/mL) | * Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges. |
|-----------|--------------------|--|
| Control 1 | 2.16-3.24 | |
| Control 2 | 51.2-76.8 | |

| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.084 |
| C | 0.252 |
| D | 0.899 |
| E | 2.82 |
| F | 9.94 |
| G | 30.0 |
| H | 76.3 |

Review/Approval

Patrizia Stadler Assoc. Director, QC *P. Stadler 30 Dec 2019*
 Name Title Signature/Date