

| | |
|-----------------|-------------|
| Product Number: | 101351 |
| Lot Number: | 502695 |
| Expiration: | 18-Jul-2021 |

| Component | Part Number | Lot Number | Concentration |
|-------------------------------|---------------|------------------|-------------------------------|
| Beads | 103611 | 031521 | 3.17X10 ⁹ beads/mL |
| Detector/Sample Diluent | 101353 | 022710 | |
| Bead Diluent | 100458 | 031128 | |
| SBG Dilution Buffer | 100376 | 029015 | |
| SBG Concentrate | 103397 | 024007 | 50nM |
| Bead Conjugation Buffer | 100453 | 028315 | |
| Bead Wash Buffer | 101352 | 028314 | |
| Bead Blocking Buffer | 100457 | 024107 | |
| Biotinylation Reaction Buffer | 100460 | 025910 | |
| EDC, 10mg | 100022 | RM202010000056 | |
| EZ-Link NHS-PEG4 Biotin | 100348 | RM2020100000049 | |
| RGP | 103159 | N/A ¹ | |
| PSA Capture Antibody | 100463 | 034006 | 5.0 mg/ml |
| PSA Detector Antibody | 100464 | 028405 | 1.1 mg/mL |
| Calibrator A-H | 100758-100765 | 027208 | |
| Control 1 | 100768 | 027209 | |
| Control 2 | 100769 | 027209 | |

¹ RGP is not Kit Lot specific.

| | Mean Range (pg/mL) | |
|-----------|--------------------|--|
| Control 1 | 2.25-3.38 | * Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges. |
| Control 2 | 47.1-70.6 | |

| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.095 |
| C | 0.281 |
| D | 0.943 |
| E | 2.95 |
| F | 9.81 |
| G | 28.5 |
| H | 73.6 |

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 15Jun2021
Signature/Date