

| | |
|-----------------|-------------|
| Product Number: | 101351 |
| Lot Number: | 502335 |
| Expiration: | 29-Aug-2020 |

| Component | Part Number | Lot Number | Concentration |
|-------------------------------|---------------|------------------|-------------------------------|
| Beads | 103611 | 005515 | 3.07X10 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 | 1.1 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) | |
|-----------|--------------------|--|
| Control 1 | 2.16-3.24 | * Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges. |
| 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

Shivani Goel
Name

QC Manager
Title

Shivani Goel 16 Apr 2020
Signature/Date