

| | |
|-----------------|-------------|
| Product Number: | 101351 |
| Lot Number: | 503251 |
| Expiration: | 15-Apr-2022 |

| Component | Part Number | Lot Number | Concentration |
|-------------------------------|---------------|------------------|---------------------------------|
| Beads | 103611 | 108830 | 2.60 X 10 ⁹ beads/mL |
| Detector/Sample Diluent | 101353 | 125711 | |
| Bead Diluent | 100458 | 122710 | |
| SBG Dilution Buffer | 100376 | 128906 | |
| SBG Concentrate | 103397 | 120035 | 50nM |
| Bead Conjugation Buffer | 100453 | 120810 | |
| Bead Wash Buffer | 101352 | 122509 | |
| Bead Blocking Buffer | 100457 | 110524 | |
| Biotinylation Reaction Buffer | 100460 | 117226 | |
| EDC, 10mg | 100022 | RM2021090000073 | |
| EZ-Link NHS-PEG4 Biotin | 100348 | RM2021070000039 | |
| RGP | 103159 | N/A ¹ | |
| PSA Capture Antibody | 100463 | 119303 | 5.1 mg/ml |
| PSA Detector Antibody | 100464 | 108106 | 1.0 mg/ml |
| Calibrator A-H | 100758-100765 | 121603 | |
| Control 1 | 100768 | 121604 | |
| Control 2 | 100769 | 121604 | |

¹ RGP is not Kit Lot specific.

| | Mean Range (pg/mL) | |
|-----------|--------------------|----------------------------------------------------------------------------------------------------------------------|
| Control 1 | 1.69-3.70 | * Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges. |
| Control 2 | 42.7-64.0 | |

Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 0.043 |
| C | 0.237 |
| D | 0.937 |
| E | 2.62 |
| F | 8.76 |
| G | 26.2 |
| H | 61.6 |

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 03 Dec 2021
Signature/Date