

| | |
|-----------------|------------|
| Product Number: | 101351 |
| Lot Number: | 501696 |
| Expiration: | 5-Nov-2019 |

| Component | Part Number | Lot Number | Concentration |
|-------------------------------|---------------|------------------|-------------------------------|
| Beads | 103206 | 823619 | 3.29X10 ⁹ beads/mL |
| Detector/Sample Diluent | 101353 | 905820 | |
| Bead Diluent | 100458 | 904217 | |
| SBG Dilution Buffer | 100376 | 902215 | |
| SBG Concentrate | 103397 | 832317 | 50nM |
| Bead Conjugation Buffer | 100453 | 830908 | |
| Bead Wash Buffer | 101352 | 903602 | |
| Bead Blocking Buffer | 100457 | 900921 | |
| Biotinylation Reaction Buffer | 100460 | 900919 | |
| EDC, 10mg | 100022 | TI276003 | |
| EZ-Link NHS-PEG4 Biotin | 100348 | TH271791 | |
| RGP | 103159 | N/A ¹ | |
| PSA Capture Antibody | 100463 | 903614 | 5.0 mg/mL |
| PSA Detector Antibody | 100464 | 831601 | 1.0 mg/mL |
| Calibrator A-H | 100758-100765 | 833801 | |
| Controls 1 and 2 | 100768-100769 | 833802 | |

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 | 2.28-3.42 | |
| Control 2 | 49.8-74.6 | |

| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.093 |
| C | 0.278 |
| D | 0.948 |
| E | 2.95 |
| F | 10.6 |
| G | 32.1 |
| H | 75.1 |

Review/Approval

Junpil Park*

QC Analyst II

 06 Mar 2019

Name

Title

Signature/Date

* On behalf of Patrizia Stadler, Assoc. Director, QC

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.