

| | |
|-----------------|-------------|
| Product Number: | 101351 |
| Lot Number: | 501297 |
| Expiration: | 17-Apr-2019 |
| Platform(s): | HD-1 |

| Component | Part Number | Lot Number | Concentration |
|-------------------------------|-----------------|------------------|-------------------------------|
| Beads | 103206 | MF-052418-01 | 3.23X10 ⁹ beads/mL |
| Detector/Sample Diluent | 101353 | MF-052318-14 | |
| Bead Diluent | 100458 | MF-052318-03 | |
| SBG Dilution Buffer | 100376 | MF-052318-12 | |
| SBG Concentrate | 100439 | 81072 | 186nM |
| Bead Conjugation Buffer | 100453 | MF-051718-04 | |
| Bead Wash Buffer | 101352 | MF-051418-09 | |
| Bead Blocking Buffer | 100457 | MF-051418-07 | |
| Biotinylation Reaction Buffer | 100460 | MF-051718-08 | |
| EDC, 10mg | 100022 | TC255240 | |
| EZ-Link NHS-PEG4 Biotin | 100348 | TC255361 | |
| RGP | 103159 | N/A ¹ | |
| PSA Capture Antibody | 100463 | 81440 | 5.1 mg/mL |
| PSA Detector Antibody | 100464 | 81441 | 1.07 mg/mL |
| Calibrator A-H | 100758-100765 | 80230 | |
| Controls 1 and 2 | 100768 & 100769 | 80231 | |

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 | 2.55-3.83 | |
| Control 2 | 57.9-88.1 | |

| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.100 |
| C | 0.327 |
| D | 1.14 |
| E | 3.44 |
| F | 12.1 |
| G | 37.1 |
| H | 93.4 |

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title


Signature/Date

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