

| | |
|------------------------|-------------|
| Product Number: | 101635 |
| Lot Number: | 501225 |
| Expiration: | 18-Oct-2018 |
| Platform(s): | HD-1 |



Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Component | Part Number | Lot Number |
|------------------------|-------------|------------------|
| Bead Reagent | 101637 | 81091 |
| Detector Reagent | 101636 | 81144 |
| SBG Reagent | 101638 | 81083 |
| Sample Diluent | 101639 | 81081 |
| Calibrator Diluent | 101640 | 81082 |
| Calibrator Concentrate | 100661 | 73463 |
| RGP Reagent | 103159 | N/A ¹ |

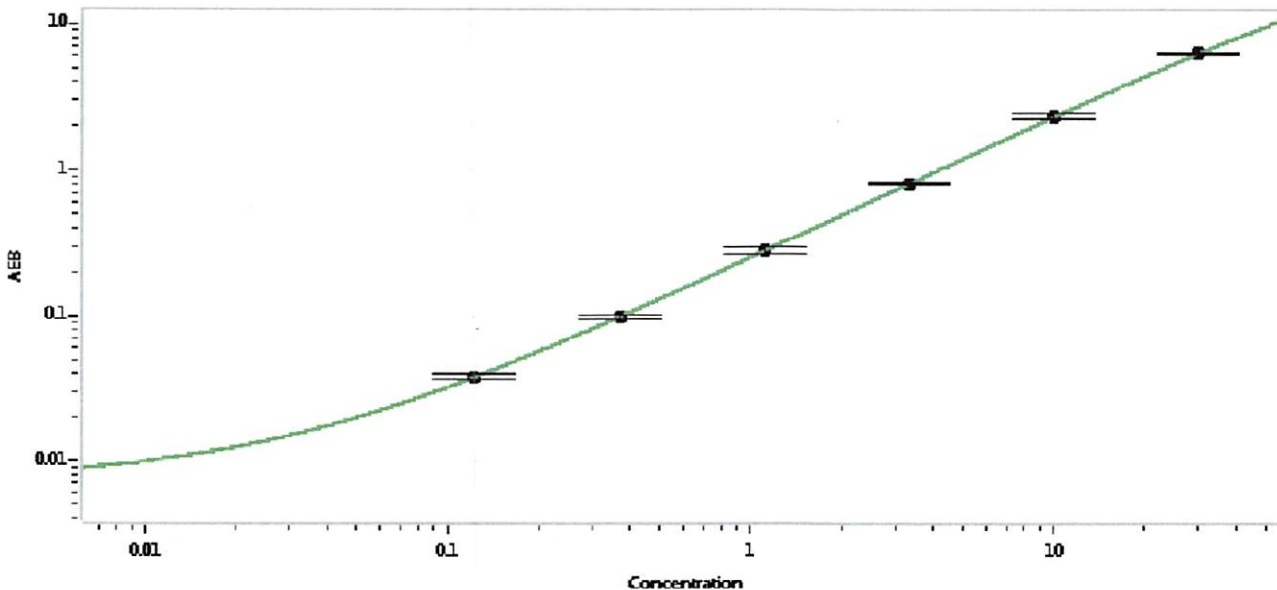
¹ RGP is not Kit Lot Specific

| Release Materials* | | | |
|--------------------|--------------|----------------|--------------------|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 52511 | 0.684 | 0.464-0.779 |
| Control 2 | 52511 | 63.4 | 43.1-78.7 |
| Panel 1 | MF-090415-01 | 0.594 | 0.355-0.780 |
| Panel 2 | MF-082515-01 | 3.73 | 2.71-4.63 |
| Panel 3 | MF-082515-02 | 35.3 | 25.0-45.3 |

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

| | |
|---|------|
| Calibrator Concentrate Value Assignment (pg/mL) | 1516 |
|---|------|

Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.123 |
| C | 0.370 |
| D | 1.11 |
| E | 3.33 |
| F | 10.0 |
| G | 30.0 |

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date
27 APR 2018