

| | |
|-----------------|-------------|
| Product Number: | 101599 |
| Lot Number: | 501355 |
| Expiration: | 26-Apr-2019 |
| Platform(s): | HD-1 & SR-X |



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

| Component | Item Number | Lot Number |
|------------------------|-------------|------------------|
| Bead Reagent | 101601 | 821302 |
| Detector Reagent | 102012 | 820704 |
| SBG Reagent | 101602 | 820705 |
| Sample Diluent | 101603 | 820708 |
| Calibrator Diluent | 102014 | 819813 |
| Calibrator Concentrate | 100778 | 81562 |
| RGP Reagent | 103159 | N/A ¹ |

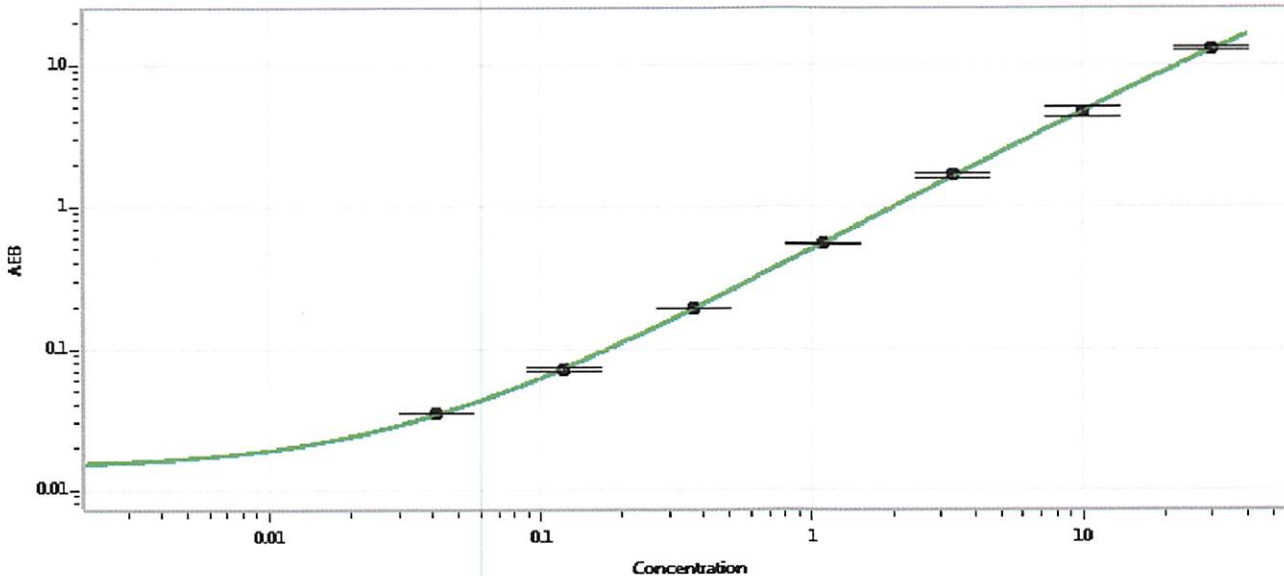
¹ RGP is not Kit Lot Specific

| Release Materials* | | | |
|--------------------|--------------|----------------|--------------------|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | MF-012518-02 | 0.464 | 0.373-0.559 |
| Control 2 | MF-012518-02 | 21.9 | 15.2-22.7 |
| Panel 1 | MF-071615-01 | 0.377 | 0.325-0.513 |
| Panel 2 | MF-071615-02 | 8.01 | 5.58-8.37 |
| Panel 3 | MF-071615-03 | 2.26 | 1.68-2.52 |

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

| | |
|---|-----|
| Calibrator Concentrate Value Assignment (pg/mL) | 800 |
|---|-----|

Calibration Curve



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 0.041 |
| C | 0.123 |
| D | 0.370 |
| E | 1.11 |
| F | 3.33 |
| G | 10.0 |
| H | 30.0 |

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date 04 Aug 2018