

|                 |            |
|-----------------|------------|
| Product Number: | 100906     |
| Lot Number:     | 501101     |
| Expiration:     | 4-Jun-2018 |
| Platform(s):    | HD-1       |



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

| Component              | Part Number | Lot Number       |
|------------------------|-------------|------------------|
| Bead Reagent           | 100911      | 73456            |
| Detector Reagent       | 100908      | 73423            |
| SBG Reagent            | 100914      | 73496            |
| Sample Diluent         | 100916      | 73388            |
| Calibrator Diluent     | 100918      | 73124            |
| Calibrator Concentrate | 100919      | 80200            |
| RGP Reagent            | 103159      | N/A <sup>1</sup> |

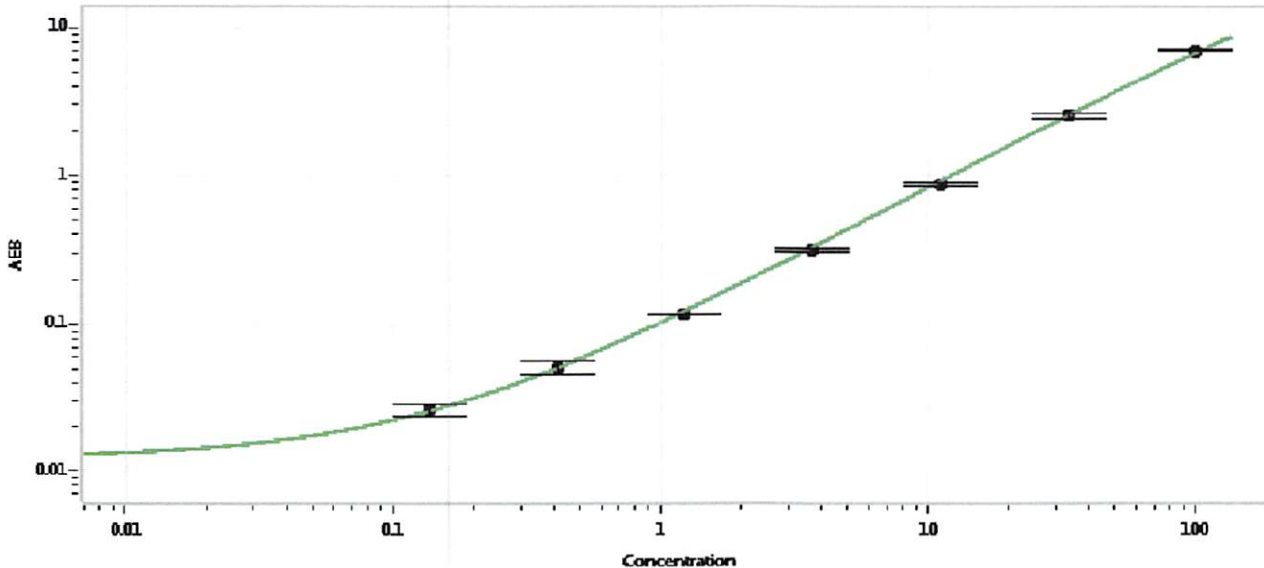
<sup>1</sup> RGP is not Kit Lot Specific

| Release Materials* |              |                |                    |
|--------------------|--------------|----------------|--------------------|
|                    | Lot          | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1          | MF-051717-02 | 1.80           | 1.52-2.54          |
| Control 2          | MF-051717-02 | 51.2           | 28.3-61.9          |
| Panel 1            | MF-032217-12 | 61.7           | 19.3-84.7          |
| Panel 2            | MF-051817-02 | 42.6           | 27.3-53.3          |
| Panel 3            | MF-051817-03 | 177            | 113-190            |

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

|   |      |
|---|------|
| Calibrator Concentrate Value Assignment (pg/mL) | 2053 |
|---|------|

Calibration Curve



Calibrator Levels (pg/mL)

|   |       |
|---|-------|
| A | 0.000 |
| B | 0.137 |
| C | 0.412 |
| D | 1.23  |
| E | 3.70  |
| F | 11.1  |
| G | 33.3  |
| H | 100   |

Review/Approval: Sivasankari Murugan (Name), QC Supervisor (Title), [Signature] (Signature), 14 FEB 2018 (Date)