

| | |
|------------------------|-----------|
| Product Number: | 102100 |
| Lot Number: | 501045 |
| Expiration: | 29-May-18 |
| Platform(s): | HD-1 |

| Component | Part Number | Lot Number |
|------------------------|-------------|--------------|
| Bead Stock | 102092 | MF-122017-02 |
| Detector Stock | 102093 | MF-011918-09 |
| Calibrator Concentrate | 102096 | MF-060817-03 |
| Calibrator Diluent | 102094 | MF-112917-17 |
| Sample Diluent | 102095 | MF-112917-19 |

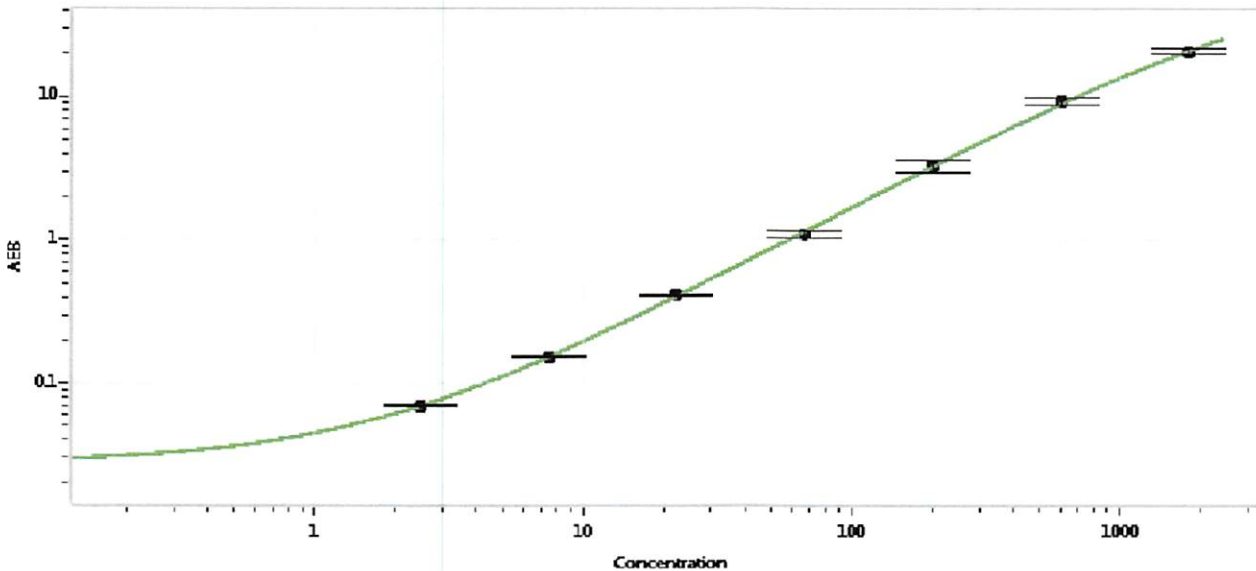


Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Generic Components - Included With Kit But Not Lot Number Specific | | |
|--|--|--------|
| Discovery Bead Diluent | | 101954 |
| Discovery Detector Diluent | | 101955 |
| Discovery SBG Diluent | | 101956 |
| Discovery SBG Concentrate | | 101957 |
| RGP Reagent | | 103159 |

| | |
|--|----------|
| Calibrator Concentrate Target Value | 36 ng/mL |
|--|----------|

Calibration Curve



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 2.47 |
| C | 7.41 |
| D | 22.2 |
| E | 66.6 |
| F | 200 |
| G | 600 |
| H | 1800 |

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 22 Jan 2018
Signature/Date