

|                 |             |
|-----------------|-------------|
| Product Number: | 101984      |
| Lot Number:     | 501471      |
| Expiration:     | 20-Feb-2019 |
| Platform(s):    | HD-1        |



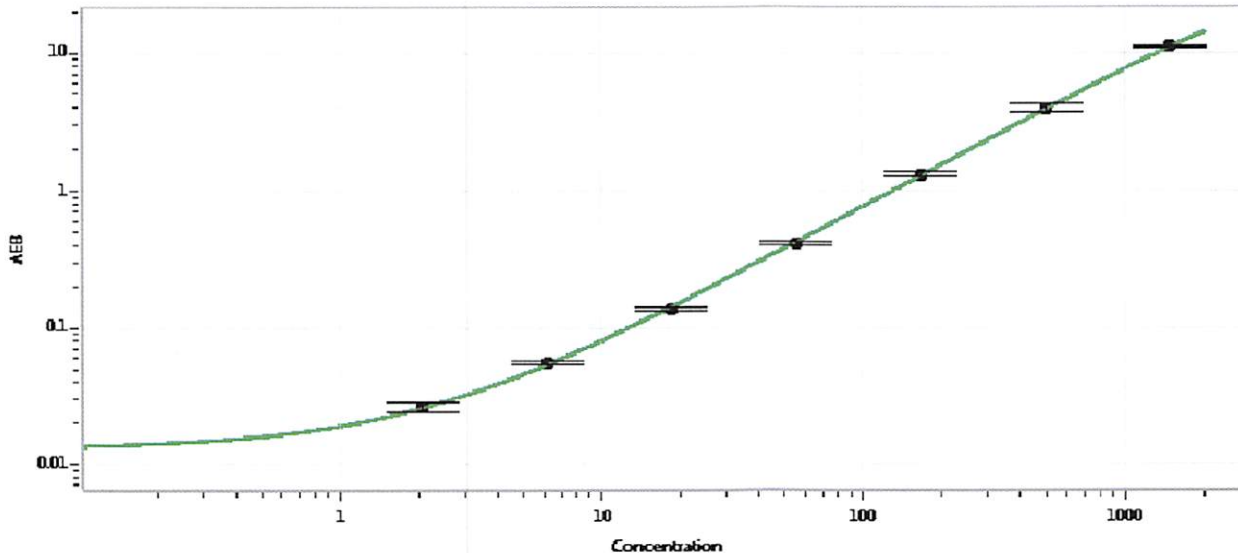
Data below represents results generated on the Simoa™ HD-1

| Component              | Part Number | Lot Number   |
|------------------------|-------------|--------------|
| Bead Stock             | 101910      | 827503       |
| Detector Stock         | 101911      | 829807       |
| Calibrator Concentrate | 101983      | MF-121417-01 |
| Calibrator Diluent     | 101908      | 826904       |
| Sample Diluent         | 101909      | 826903       |

| Generic Components - Included With Kit But Not Lot Number Specific |        |
|--|--------|
| Discovery Bead Diluent   | 101954 |
| Discovery Detector Diluent   | 101955 |
| Discovery SBG Diluent  | 101956 |
| Discovery SBG Stock  | 103297 |
| RGP Reagent  | 103159 |

|  |          |
|--|----------|
| Calibrator Concentrate<br>Target Value | 30 ng/mL |
|--|----------|

Calibration Curve



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 2.06  |
| C                         | 6.17  |
| D                         | 18.5  |
| E                         | 55.6  |
| F                         | 167   |
| G                         | 500   |
| H                         | 1500  |

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 06 Nov 2018  
Signature/Date