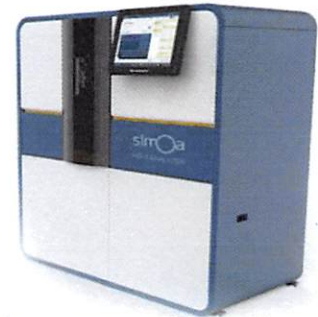


|                 |            |
|-----------------|------------|
| Product Number: | 102336     |
| Lot Number:     | 501277     |
| Expiration:     | 9-Nov-2018 |
| Platform(s):    | HD-1       |



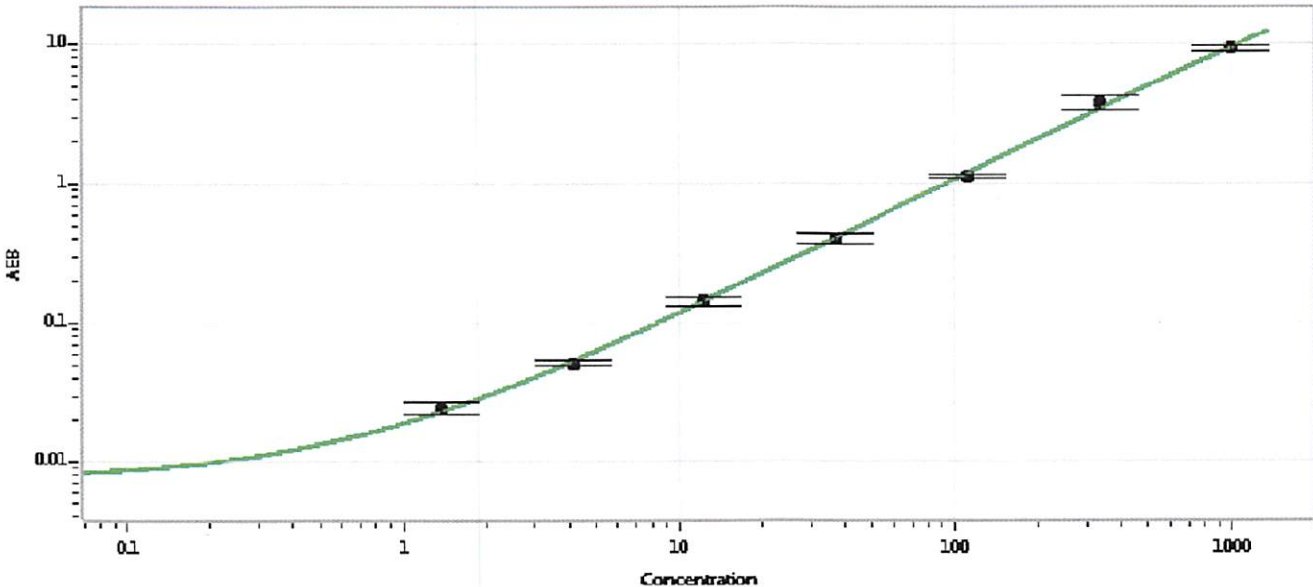
Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Component              | Part Number | Lot Number   |
|------------------------|-------------|--------------|
| Bead Stock             | 102330      | MF-052418-02 |
| Detector Stock         | 102331      | MF-052318-08 |
| Calibrator Concentrate | 102335      | MF-092917-03 |
| Calibrator Diluent     | 102332      | MF-050918-16 |
| Sample Diluent         | 102333      | MF-052318-09 |

| 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 | 20 ng/mL |
|-------------------------------------|----------|

Calibration Curve



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 1.37  |
| C                         | 4.11  |
| D                         | 12.3  |
| E                         | 37.0  |
| F                         | 111   |
| G                         | 333   |
| H                         | 1000  |

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date  
08 Nov 2018