

|                 |             |
|-----------------|-------------|
| Product Number: | 103093      |
| Lot Number:     | 501356      |
| Expiration:     | 15-Oct-2018 |
| Platform(s):    | HD-1 & SR-X |

| Component              | Part Number | Lot Number   |
|------------------------|-------------|--------------|
| Bead Stock             | 102021      | MF-041518-02 |
| Detector Stock         | 103043      | 821310       |
| Calibrator Concentrate | 103047      | MF-010818-10 |
| Calibrator Diluent     | 103044      | MF-041918-04 |
| Sample Diluent         | 103046      | MF-041918-05 |

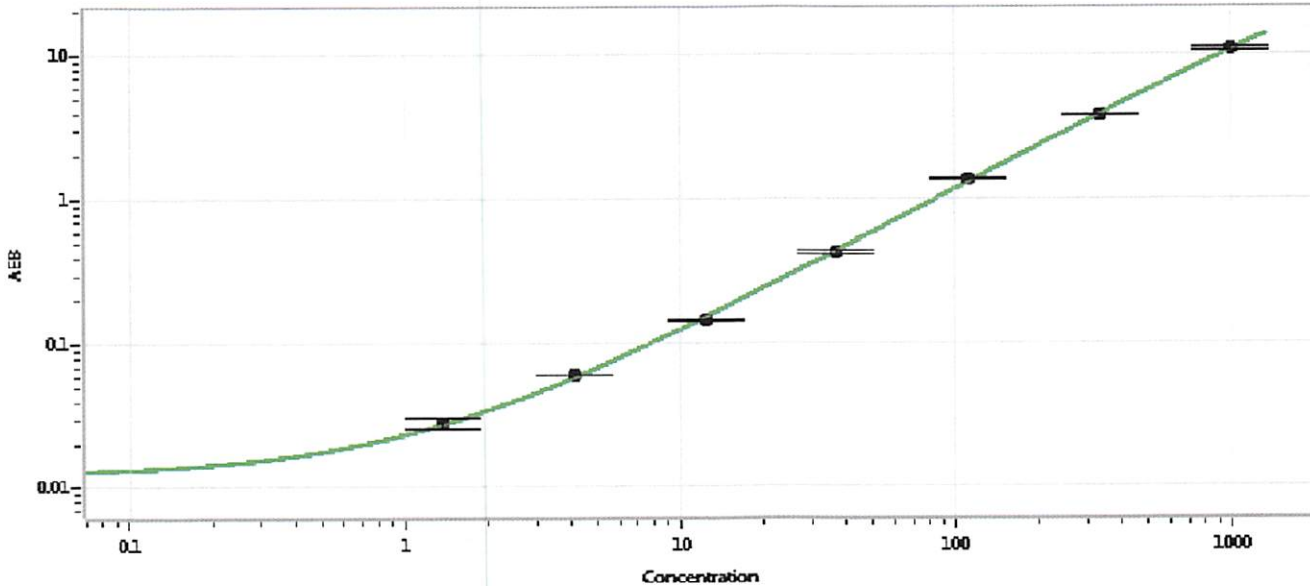


Data below represents results generated on the Simoa™ HD-1 Analyzer.

| 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 | 12 ng/mL |
|-------------------------------------|----------|

Calibration Curve



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 1.37  |
| C                         | 4.12  |
| D                         | 12.3  |
| E                         | 37.0  |
| F                         | 111   |
| G                         | 333   |
| H                         | 1000  |

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date  
02 AUG 2018