

|                 |             |
|-----------------|-------------|
| Product Number: | 102336      |
| Lot Number:     | 501950      |
| Expiration:     | 30-Mar-2020 |
| Platform(s):    | HD-1 / HD-X |



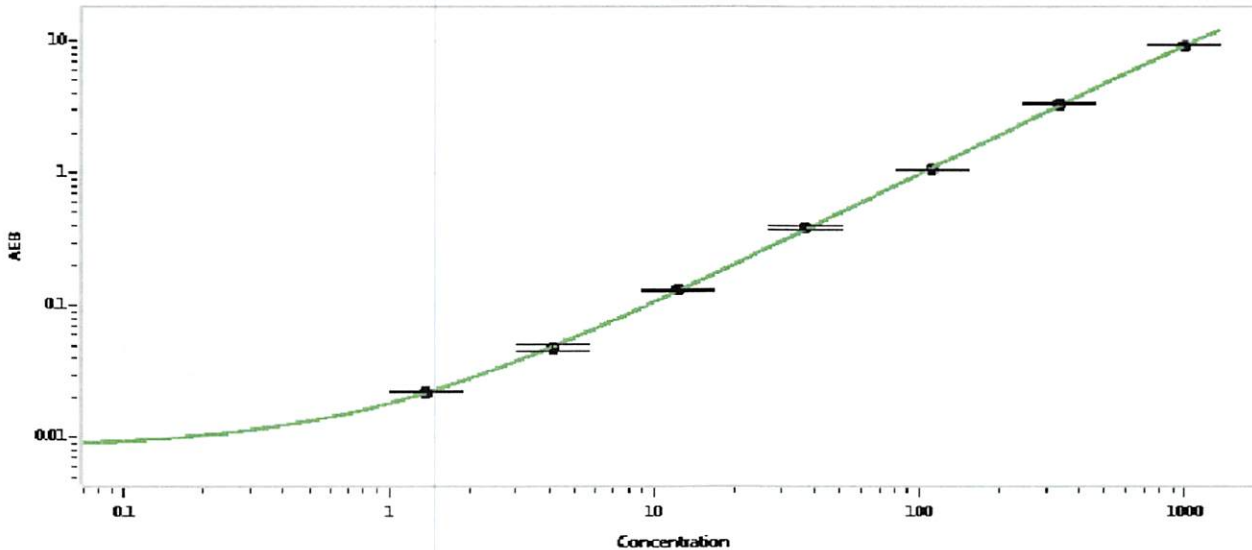
Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Component              | Part Number | Lot Number |
|------------------------|-------------|------------|
| Bead Stock             | 102330      | 920309     |
| Detector Stock         | 102331      | 920303     |
| Calibrator Concentrate | 102335      | 920503     |
| Calibrator Diluent     | 102332      | 908917     |
| Sample Diluent         | 102333      | 920304     |


| 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 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 | *Junpil Park  | QC Analyst II |  | 30 Jul 2019 |
|                 | Name  | Title         | Signature/Date  |             |
|                 | *On behalf of Patrizia Stadler, Assoc. Director, QC |               |   |             |