

|                 |             |
|-----------------|-------------|
| Product Number: | 102635      |
| Lot Number:     | 501884      |
| Expiration:     | 27-Feb-2020 |
| Platform(s):    | HD-1/ HD-X  |



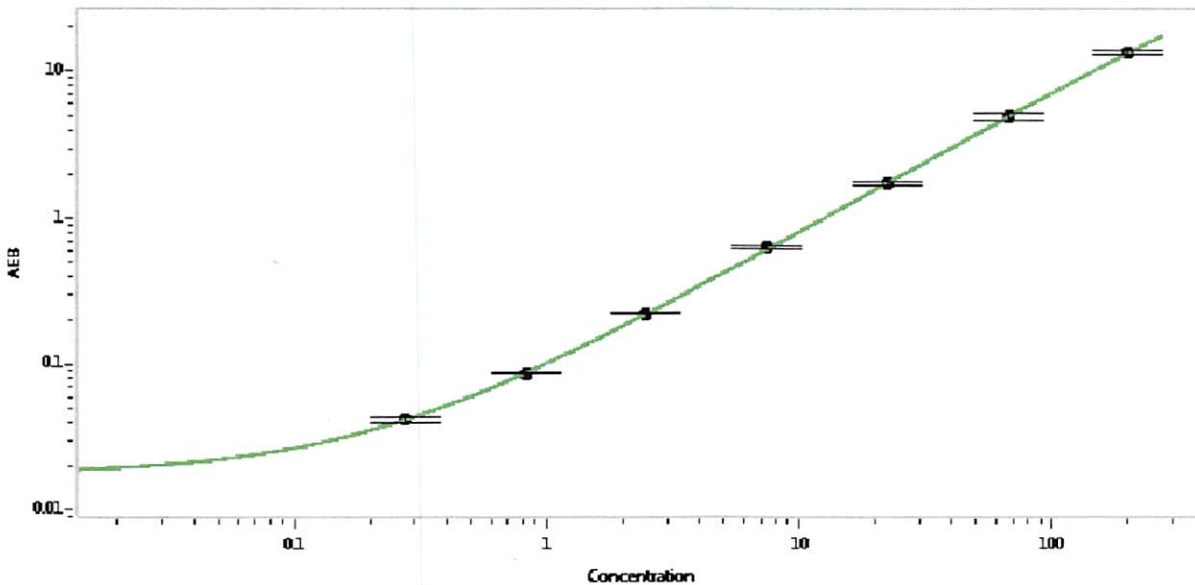
Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Component              | Part Number | Lot Number |
|------------------------|-------------|------------|
| Bead Stock             | 102628      | 909210     |
| Detector Stock         | 102629      | 909309     |
| Calibrator Concentrate | 102632      | 822101     |
| Calibrator Diluent     | 102630      | 905825     |
| Sample Diluent         | 102631      | 909310     |

| 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 | 4 ng/mL |
|-------------------------------------|---------|

Calibration Curve



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 0.274 |
| C                         | 0.823 |
| D                         | 2.47  |
| E                         | 7.41  |
| F                         | 22.2  |
| G                         | 66.7  |
| H                         | 200   |

|                 |                  |                     |                               |
|-----------------|------------------|---------------------|-------------------------------|
| Review/Approval | Patrizia Stadler | Assoc. Director, QC | <i>P. Stadler</i> 27 JUN 2019 |
|                 | Name             | Title               | Signature/Date                |