

| | |
|-----------------|-------------|
| Product Number: | 101871 |
| Lot Number: | 502455 |
| Expiration: | 22-Nov-2020 |
| Platform(s): | HD-1/HD-X |



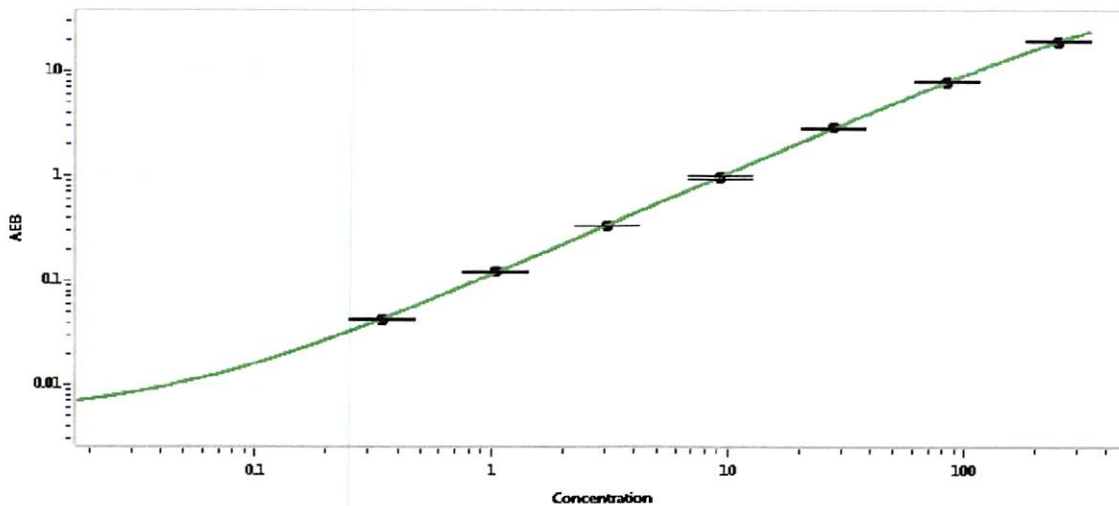
Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Component | Part Number | Lot Number |
|------------------------|-------------|------------|
| Bead Stock | 101893 | 001605 |
| Detector Stock | 101894 | 001601 |
| Calibrator Concentrate | 101866 | 926009 |
| Calibrator Diluent | 101959 | 001316 |
| Sample Diluent | 101958 | 932625 |

| 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 | 5 ng/mL |
|-------------------------------------|---------|

Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.343 |
| C | 1.03 |
| D | 3.09 |
| E | 9.26 |
| F | 27.8 |
| G | 83.3 |
| H | 250 |

Review/Approval

Shivani Goel

Quality Control Manager

Shivani Goel 06/11/2020

Name

Title

Signature/Date