

| | |
|-----------------|-------------|
| Product Number: | 102336 |
| Lot Number: | 502185 |
| Expiration: | 21-Nov-2020 |
| Platform(s): | HD-1/HD-X |



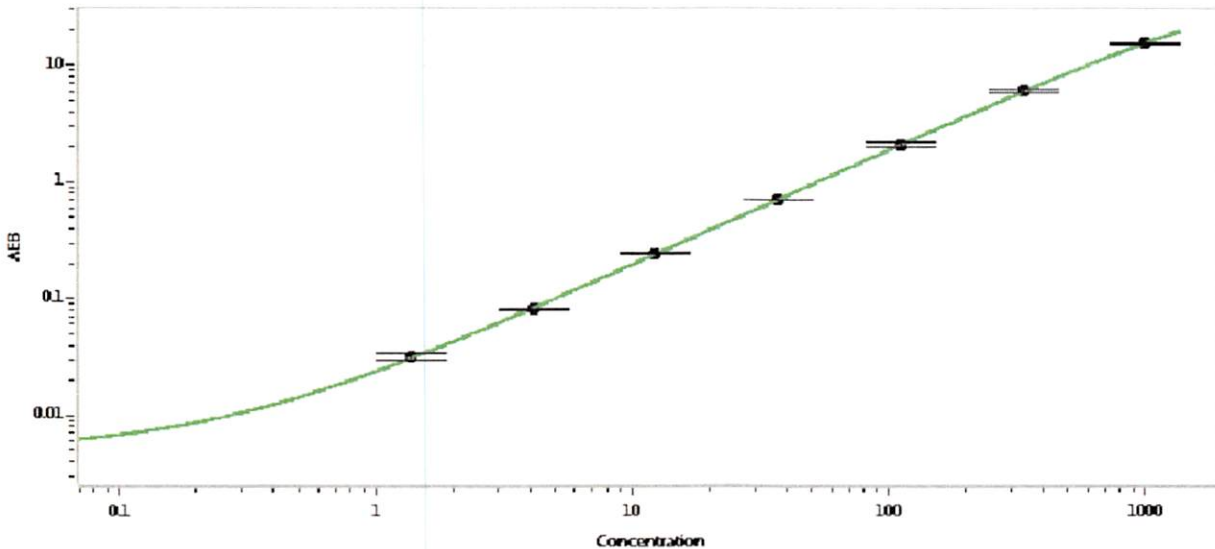
Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Component | Part Number | Lot Number |
|------------------------|-------------|------------|
| Bead Stock | 102330 | 932914 |
| Detector Stock | 102331 | 932613 |
| Calibrator Concentrate | 102335 | 933104 |
| Calibrator Diluent | 102332 | 932615 |
| Sample Diluent | 102333 | 932617 |

| 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

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 09 Jan 2020
Signature/Date