

| | |
|-----------------|-------------|
| Product Number: | 101984 |
| Lot Number: | 502074 |
| Expiration: | 29-Jul-2020 |
| Platform(s): | HD-1/HD-X |

| Component | Part Number | Lot Number |
|------------------------|-------------|------------|
| Bead Stock | 101910 | 928103 |
| Detector Stock | 101911 | 928010 |
| Calibrator Concentrate | 101983 | 910010 |
| Calibrator Diluent | 101908 | 928203 |
| Sample Diluent | 101909 | 928011 |

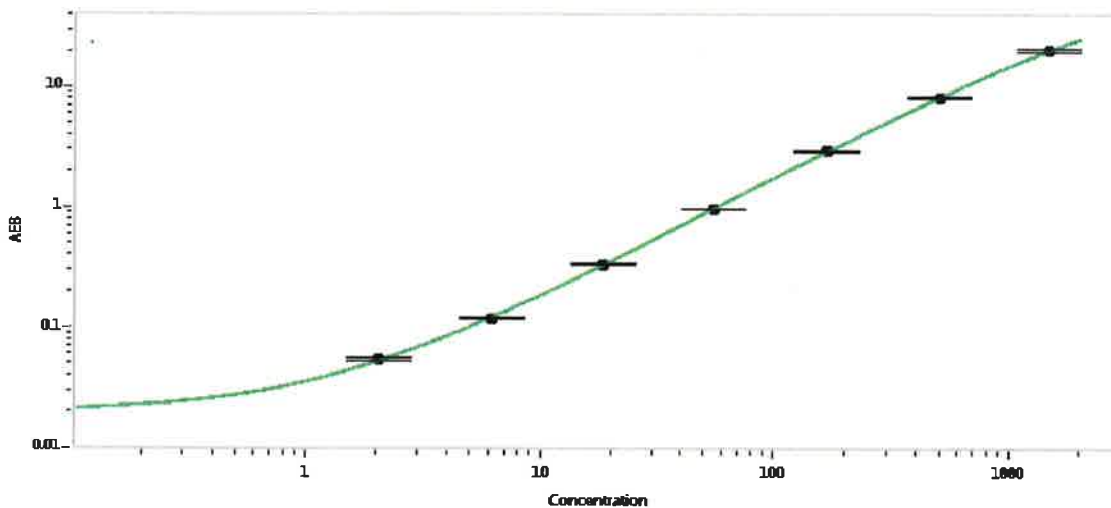


Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Generic Components - Included With Kit But Not Lot Number Specific | |
|--|--------|
| Discovery Bead Diluent | 101954 |
| Discovery Detector Diluent | 101955 |
| Discovery SBG Diluent | 101956 |
| Discovery SBG Stock | 103297 |
| RGP Reagent | 103159 |

| | |
|-------------------------------------|----------|
| Calibrator Concentrate Target Value | 30 ng/mL |
|-------------------------------------|----------|

Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 2.06 |
| C | 6.17 |
| D | 18.5 |
| E | 55.6 |
| F | 167 |
| G | 500 |
| H | 1500 |

| | | | |
|-----------------|------------------|---------------------|-------------------------------|
| Review/Approval | Patrizia Stadler | Assoc. Director, QC | <i>P. Stadler</i> 23 Oct 2019 |
| | Name | Title | Signature/Date |