

| | |
|-----------------|------------------|
| Product Number: | 102209 |
| Lot Number: | 501969 |
| Expiration: | 19-Jun-2020 |
| Platform(s): | HD-1/HD-X & SR-X |



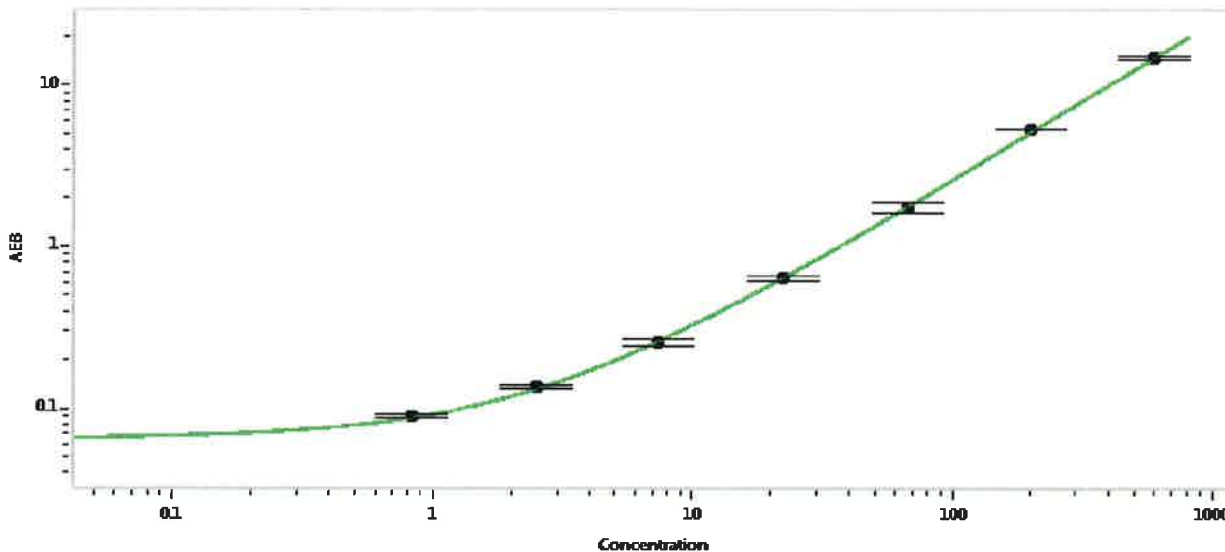
Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Component | Part Number | Lot Number |
|------------------------|-------------|------------|
| Bead Stock | 102201 | 921405 |
| Detector Stock | 102202 | 921309 |
| Calibrator Concentrate | 102205 | 827108 |
| Calibrator Diluent | 102203 | 921207 |
| Sample Diluent | 102204 | 921208 |

| 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 | 12 ng/mL |
|-------------------------------------|----------|

Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.823 |
| C | 2.47 |
| D | 7.41 |
| E | 22.2 |
| F | 66.7 |
| G | 200 |
| H | 600 |

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 27 Aug 2019
Signature/Date