

| | |
|-----------------|-------------|
| Product Number: | 100860 |
| Lot Number: | 501928 |
| Expiration: | 23-Apr-2020 |
| Platform(s): | HD-1/ HD-X |



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

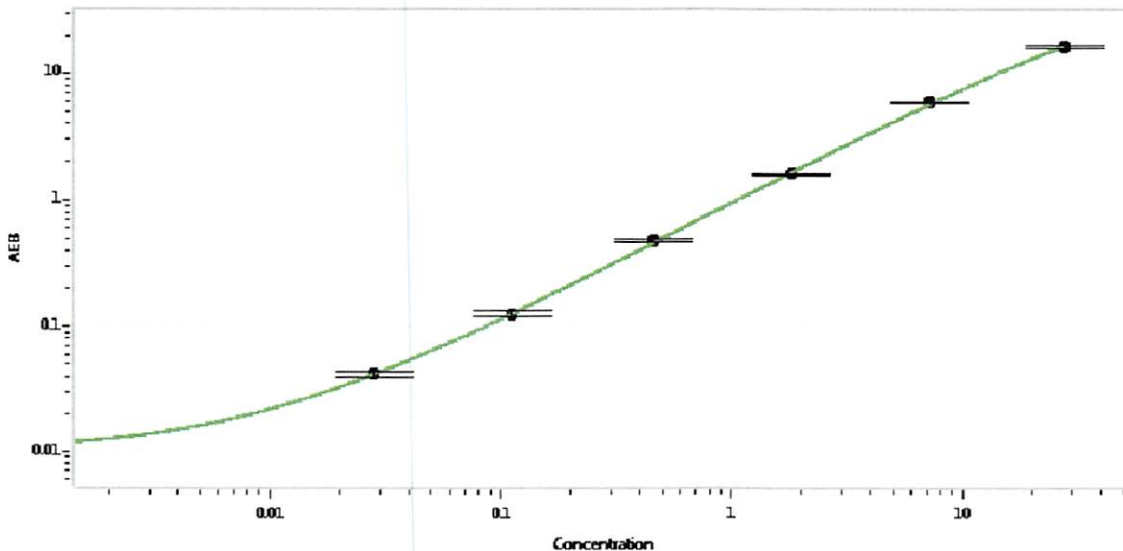
| Component | Part Number | Lot Number |
|------------------|---------------|------------------|
| Bead Reagent | 100865 | 911410 |
| Detector Reagent | 100862 | 911506 |
| SBG Reagent | 100868 | 911311 |
| Sample Diluent | 100870 | 911310 |
| Calibrators | 101111 | 907103 |
| Controls | 100875-100876 | 910602 |
| RGP Reagent | 103159 | N/A ¹ |

¹ RGP is not Kit Lot Specific

| Release Materials* | | | |
|--------------------|--------|----------------|--------------------|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 910602 | 6.11 | 3.46-10.4 |
| Control 2 | 910602 | 0.110 | 0.060-0.180 |
| Panel 1 | 908501 | 1.40 | 1.27-1.91 |
| Panel 2 | 908502 | 9.65 | 7.66-12.0 |
| Panel 3 | 908503 | 34.1 | 24.2-45.5 |

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.028 |
| C | 0.112 |
| D | 0.459 |
| E | 1.82 |
| F | 7.20 |
| G | 27.30 |

| | | | |
|-----------------|------------------|---------------------|-------------------------------|
| Review/Approval | Patrizia Stadler | Assoc. Director, QC | <i>P. Stadler</i> 25 Jul 2019 |
| | Name | Title | Signature/Date |