

| | |
|-----------------|------------|
| Product Number: | 100860 |
| Lot Number: | 501744 |
| Expiration: | 6-Mar-2020 |
| Platform(s): | HD-1 |



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

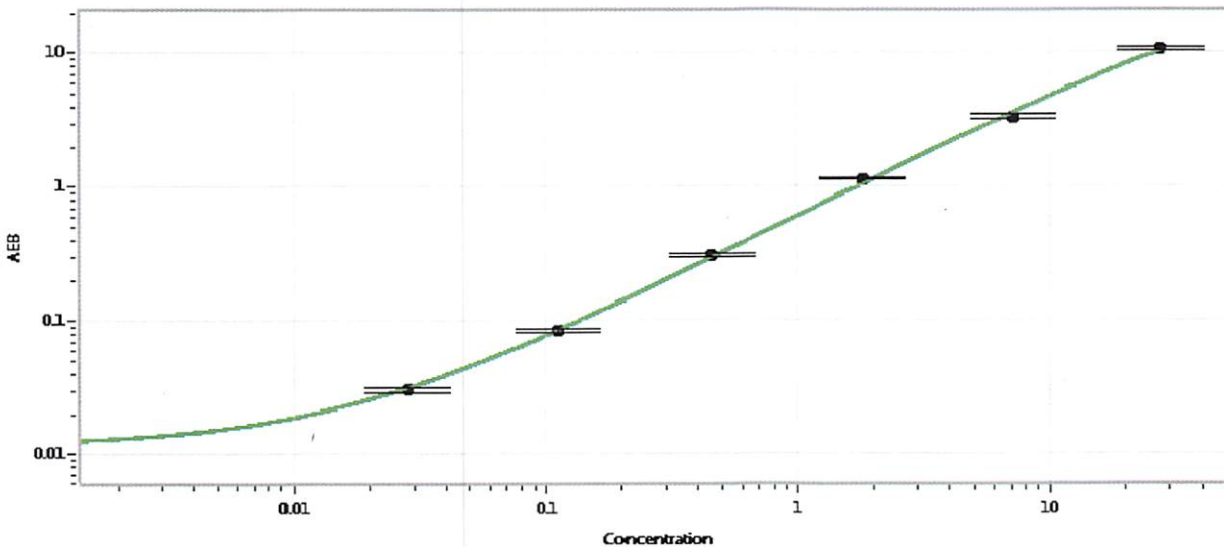
| Component | Part Number | Lot Number |
|------------------|-----------------|------------------|
| Bead Reagent | 100865 | 908706 |
| Detector Reagent | 100862 | 906508 |
| SBG Reagent | 100868 | 906507 |
| Sample Diluent | 100870 | 906506 |
| Calibrators | 101111 | 907103 |
| Controls | 100875 & 100876 | 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 | 6.29 | 3.53-10.6 |
| Control 2 | 910602 | 0.117 | 0.061-0.182 |
| Panel 1 | 908501 | 1.68 | 1.30-1.94 |
| Panel 2 | 908502 | 9.37 | 7.54-12.2 |
| Panel 3 | 908503 | 39.0 | 23.6-46.4 |

* 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 | 27.3 |

| | | | | |
|-----------------|------------------|---------------------------|-------------------|-------------|
| Review/Approval | Patrizia Stadler | Associate Director, QC | <i>P. Stodler</i> | 22 Apr 2019 |
| | Name | Title | Signature/Date | |