

| | |
|-----------------|-------------|
| Product Number: | 100196 |
| Lot Number: | 502418 |
| Expiration: | 12-Mar-2021 |
| Platform(s): | HD-1/HD-X |



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

| Component | Part Number | Lot Number |
|------------------------|-------------|------------------|
| Bead Reagent | 100391 | 016306 |
| Detector Reagent | 100315 | 016109 |
| SBG Reagent | 100316 | 016002 |
| Sample Diluent | 100317 | 015414 |
| Calibrator Diluent | 100665 | 016102 |
| Calibrator Concentrate | 100664 | 907102 |
| RGP Reagent | 103159 | N/A ¹ |

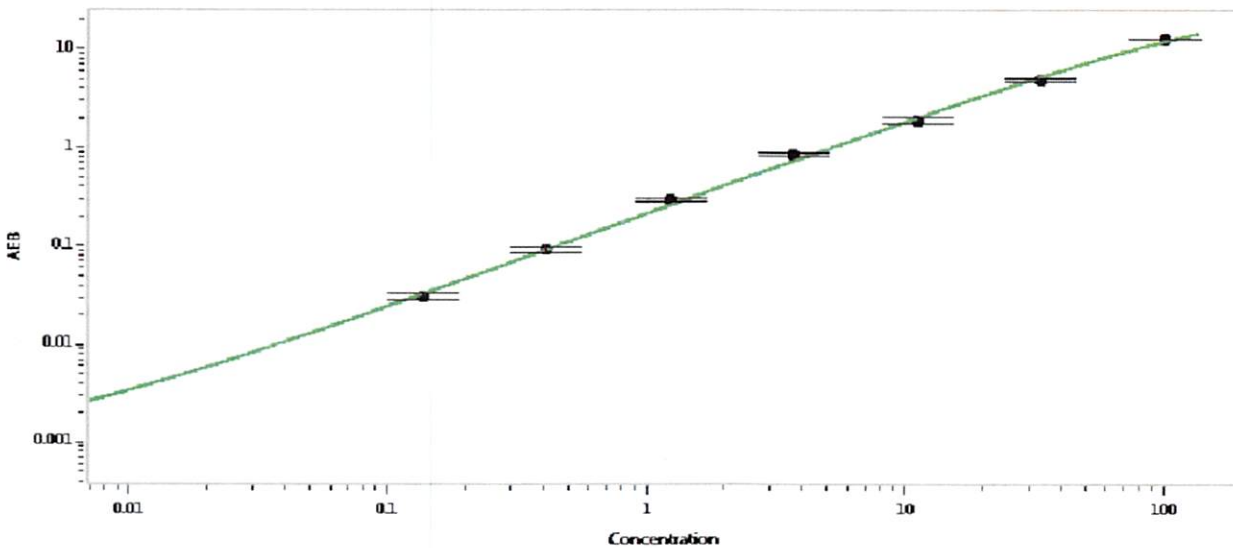
¹ RGP is not Kit Lot Specific

| Release Materials* | | | |
|--------------------|--------------|----------------|--------------------|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 910002 | 9.50 | 7.24-10.9 |
| Control 2 | 910002 | 86.9 | 73.9-111 |
| Panel 1a | MF-062018-01 | 0.314 | 0.097-1.20 |
| Panel 2 | MF-062018-02 | 8.74 | 6.90-10.9 |
| Panel 3 | MF-062018-03 | 91.0 | 72.9-155 |

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

| | |
|---|------|
| Calibrator Concentrate Value Assignment (pg/mL) | 2290 |
|---|------|

Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.137 |
| C | 0.412 |
| D | 1.23 |
| E | 3.70 |
| F | 11.1 |
| G | 33.3 |
| H | 100 |

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 02/11/2020
Signature/Date