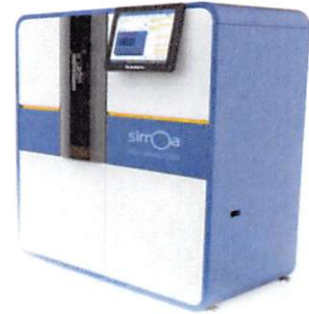


|                        |             |
|------------------------|-------------|
| <b>Product Number:</b> | 101635      |
| <b>Lot Number:</b>     | 501805      |
| <b>Expiration:</b>     | 29-Mar-2020 |
| <b>Platform(s):</b>    | HD-1        |



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

| Component              | Part Number | Lot Number       |
|------------------------|-------------|------------------|
| Bead Reagent           | 101637      | 910811           |
| Detector Reagent       | 101636      | 909506           |
| SBG Reagent            | 101638      | 912012           |
| Sample Diluent         | 101639      | 908813           |
| Calibrator Diluent     | 101640      | 908812           |
| Calibrator Concentrate | 100661      | 907709           |
| RGP Reagent            | 103159      | N/A <sup>1</sup> |

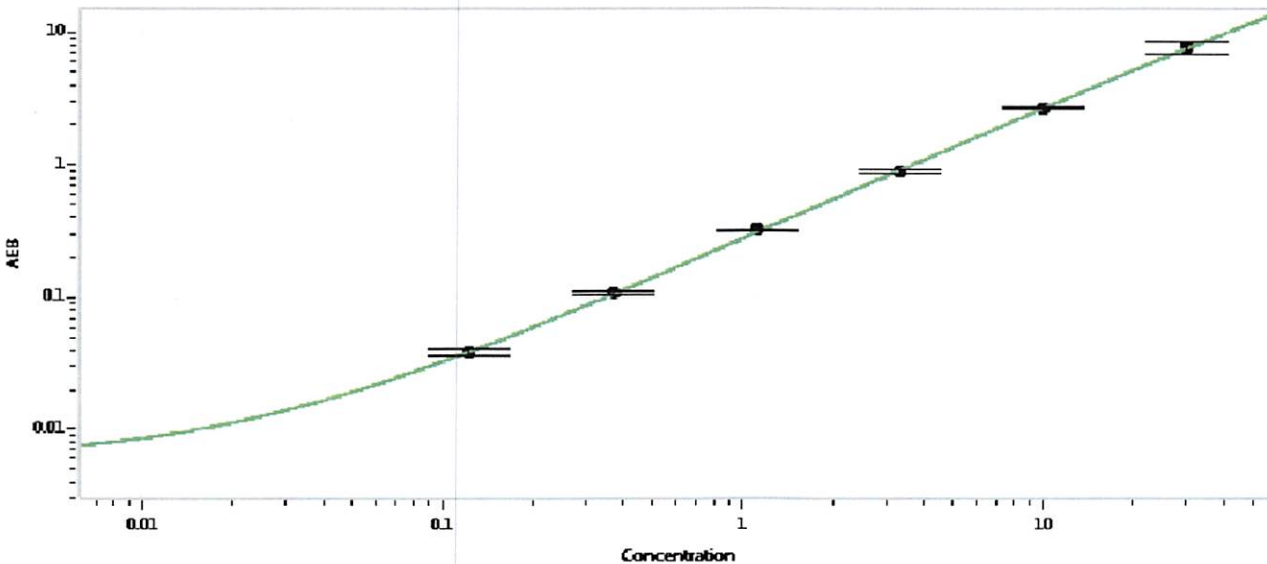
<sup>1</sup> RGP is not Kit Lot Specific

| Release Materials* |              |                |                    |
|--------------------|--------------|----------------|--------------------|
|                    | Lot          | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1          | 904406       | 0.734          | 0.558-0.837        |
| Control 2          | 904406       | 71.6           | 53.6-80.4          |
| Panel 1            | MF-040618-02 | 0.540          | 0.407-0.654        |
| Panel 2            | MF-040618-03 | 5.66           | 4.18-6.27          |
| Panel 3            | MF-040618-04 | 54.0           | 43.8-65.7          |

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

|   |      |
|---|------|
| Calibrator Concentrate Value Assignment (pg/mL) | 1810 |
|---|------|

Calibration Curve



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 0.123 |
| C                         | 0.370 |
| D                         | 1.11  |
| E                         | 3.33  |
| F                         | 10.0  |
| G                         | 30.0  |

Review/Approval: Patrizia Stadler Assoc. Director, QC  
 Name: Patrizia Stadler Title: Assoc. Director, QC  
 Signature/Date: P. Stadler 06 MAY 2019