

|                 |            |
|-----------------|------------|
| Product Number: | 100196     |
| Lot Number:     | 503733     |
| Expiration:     | 6-Aug-2023 |
| 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      | 230109           |
| Detector Reagent       | 100315      | 230405           |
| SBG Reagent            | 100316      | 230501           |
| Sample Diluent         | 100317      | 230113           |
| Calibrator Diluent     | 100665      | 230620           |
| Calibrator Concentrate | 100664      | 121803           |
| RGP Reagent            | 103159      | N/A <sup>1</sup> |

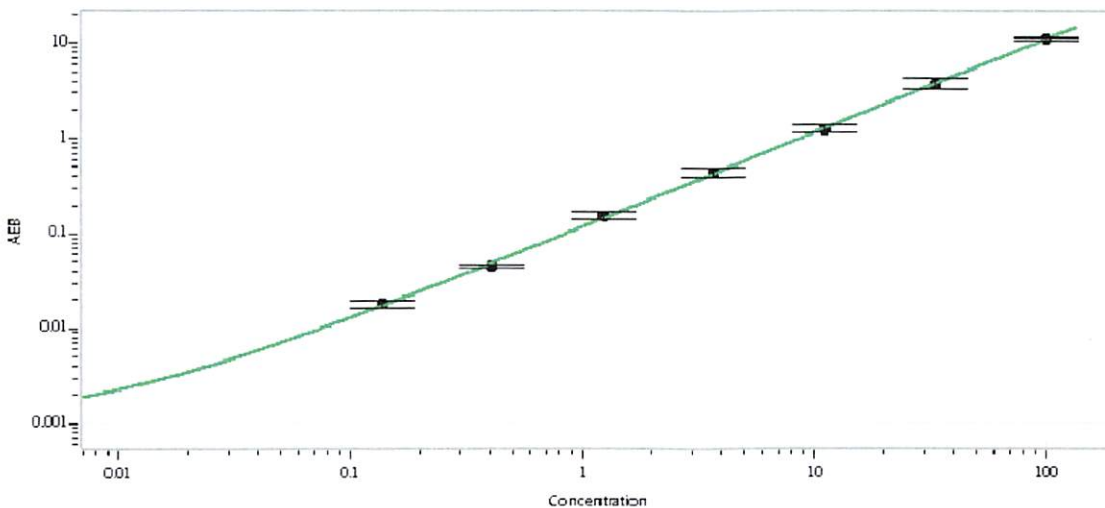
<sup>1</sup> RGP is not Kit Lot Specific

| Release Materials* |        |                |                     |
|--------------------|--------|----------------|---------------------|
|                    | Lot    | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1          | 123410 | 12.2           | 5.34 - 13.4         |
| Control 2          | 123410 | 112            | 51.9 - 143          |
| Panel 1            | 029002 | 2.43           | 1.19 - 2.82         |
| Panel 2            | 029002 | 12.1           | 5.13 - 14.0         |
| Panel 3            | 029002 | 92.2           | 46.8 - 134          |

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

|   |      |
|---|------|
| Calibrator Concentrate Value Assignment (pg/mL) | 2110 |
|---|------|

Calibration Curve



Calibrator Levels (pg/mL)

|   |       |
|---|-------|
| A | 0.000 |
| B | 0.137 |
| C | 0.412 |
| D | 1.24  |
| E | 3.70  |
| F | 11.1  |
| G | 33.3  |
| H | 100   |

Review/Approval

Tony Lem  
Name

QC  
Title

*[Signature]* 30NOV2022  
Signature/Date