

|                 |                  |
|-----------------|------------------|
| Product Number: | 103186           |
| Lot Number:     | 502034           |
| Expiration:     | 15-Apr-2020      |
| Platform(s):    | HD-1/HD-x & SR-X |



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

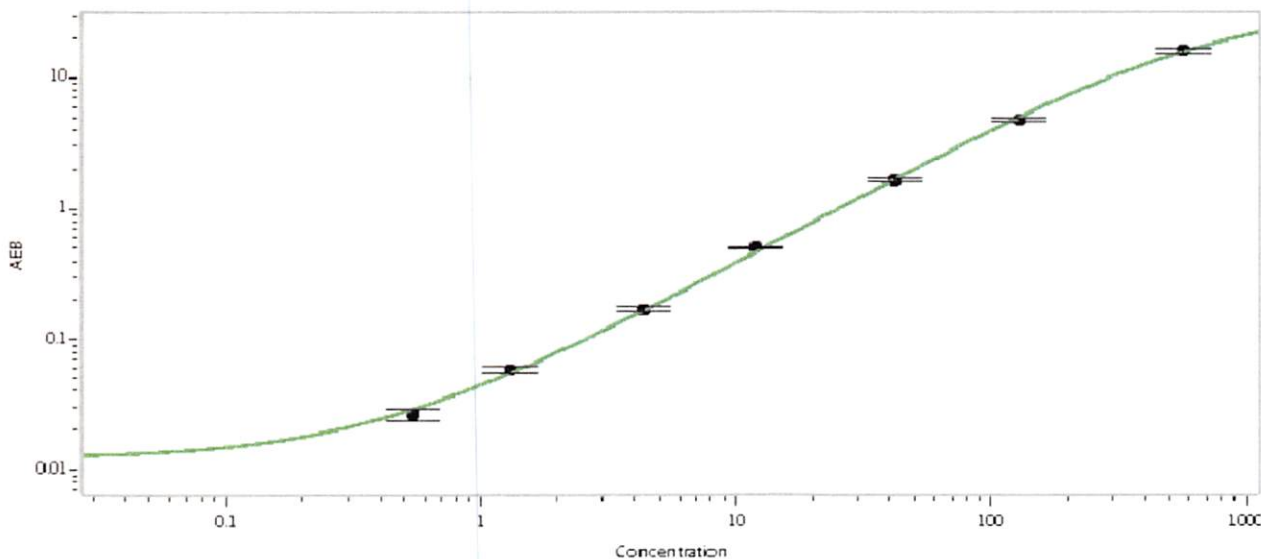
| Component        | Part Number     | Lot Number       |
|------------------|-----------------|------------------|
| Bead Reagent     | 102246          | 911413           |
| Detector Reagent | 102248          | 910508           |
| SBG Reagent      | 102250          | 910509           |
| Sample Diluent   | 102252          | 913002           |
| Calibrators      | 102255          | 925601           |
| Controls         | 102256 - 102257 | 925901           |
| 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          | 924803 | 4.09           | 3.04-4.63          |
| Control 2          | 924803 | 144            | 110-165            |
| Panel 1a           | 823301 | 5.61           | 4.08-6.68          |
| Panel 1b           | 823302 | 6.41           | 4.90-7.61          |
| Panel 2            | 823601 | 78.5           | 59.3-101           |
| Panel 3            | 823602 | 286            | 236-402            |

\* 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                         | 0.542 |
| C                         | 1.31  |
| D                         | 4.34  |
| E                         | 12.0  |
| F                         | 41.8  |
| G                         | 129   |
| H                         | 567   |

|                 |                  |                     |                    |
|-----------------|------------------|---------------------|--------------------|
| Review/Approval | Patrizia Stadler | Assoc. Director, QC | <i>p. Stadler</i>  |
|                 | Name             | Title               | Signature/Date     |
|                 |                  |                     | <i>23 Sep 2019</i> |