

|                 |                  |
|-----------------|------------------|
| Product Number: | 103186           |
| Lot Number:     | 502035           |
| Expiration:     | 16-Feb-2020      |
| Platform(s):    | HD-1/HD-x & SR-X |



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

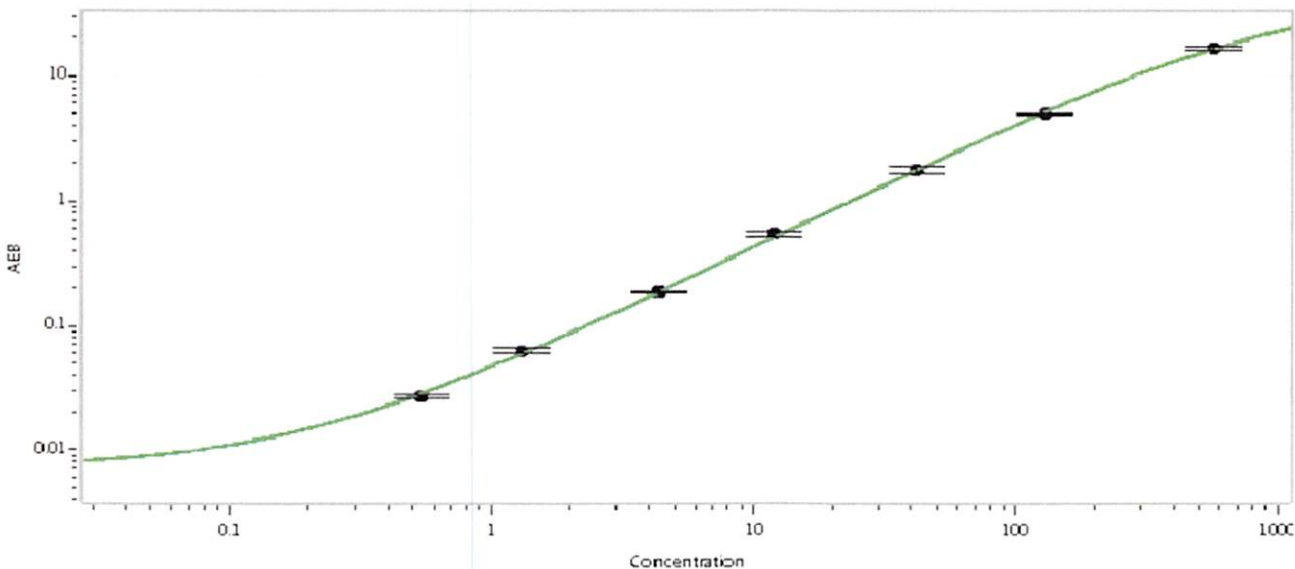
| Component        | Part Number     | Lot Number       |
|------------------|-----------------|------------------|
| Bead Reagent     | 102246          | 906503           |
| Detector Reagent | 102248          | 904703           |
| SBG Reagent      | 102250          | 904702           |
| Sample Diluent   | 102252          | 904701           |
| 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 | 3.93           | 3.04-4.63          |
| Control 2          | 924803 | 147            | 110-165            |
| Panel 1a           | 823301 | 5.52           | 4.08-6.68          |
| Panel 1b           | 823302 | 6.29           | 4.90-7.61          |
| Panel 2            | 823601 | 76.8           | 59.3-101           |
| Panel 3            | 823602 | 284            | 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 *P. Stadler* 23 Sep 2019  
Name Title Signature/Date