

|                 |             |
|-----------------|-------------|
| Product Number: | 103186      |
| Lot Number:     | 503469      |
| Expiration:     | 3-May-2023  |
| Platform(s):    | HD-1 / HD-X |



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

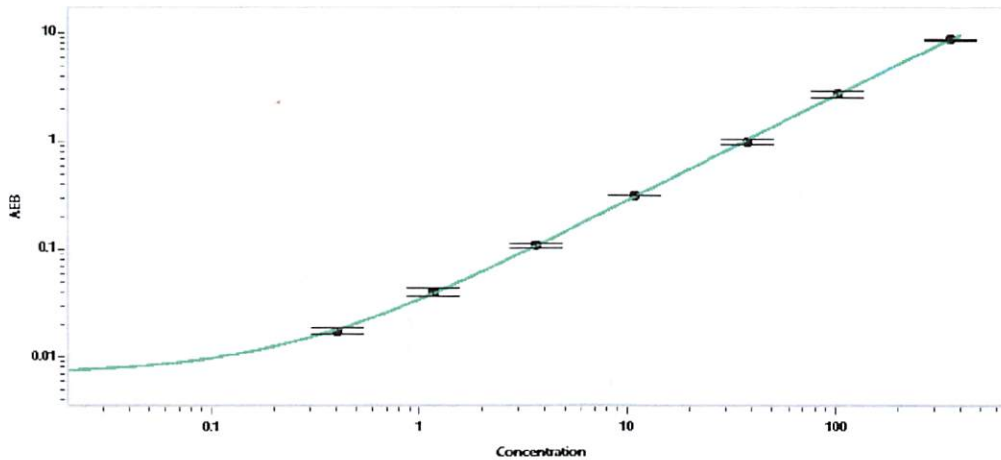
| Component        | Part Number | Lot Number       |
|------------------|-------------|------------------|
| Bead Reagent     | 102246      | 214407           |
| Detector Reagent | 102248      | 214201           |
| SBG Reagent      | 102250      | 214202           |
| Sample Diluent   | 102252      | 214203           |
| Calibrators      | 102255      | 213603           |
| Control 1        | 102256      | 213604           |
| Control 2        | 102257      | 213604           |
| 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          | 213604 | 3.80           | 2.17 - 4.36        |
| Control 2          | 213604 | 142            | 103 - 155          |
| Panel 1            | 134208 | 7.89           | 3.95 - 9.43        |
| Panel 2            | 134208 | 5.85           | 1.86 - 7.15        |
| Panel 3            | 212511 | 85.5           | 73.5 - 116         |
| Panel 4            | 212511 | 220            | 145 - 372          |

\* 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                         | 0.407 |
| C                         | 1.17  |
| D                         | 3.62  |
| E                         | 10.8  |
| F                         | 37.7  |
| G                         | 103   |
| H                         | 359   |

Review/Approval           Matt Smart          Director, QC                     [Signature]           15 Jun 2022  
 Name Title Signature/Date