

|                 |            |
|-----------------|------------|
| Product Number: | 100860     |
| Lot Number:     | 501595     |
| Expiration:     | 9-Jul-2019 |
| Platform(s):    | HD-1       |



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

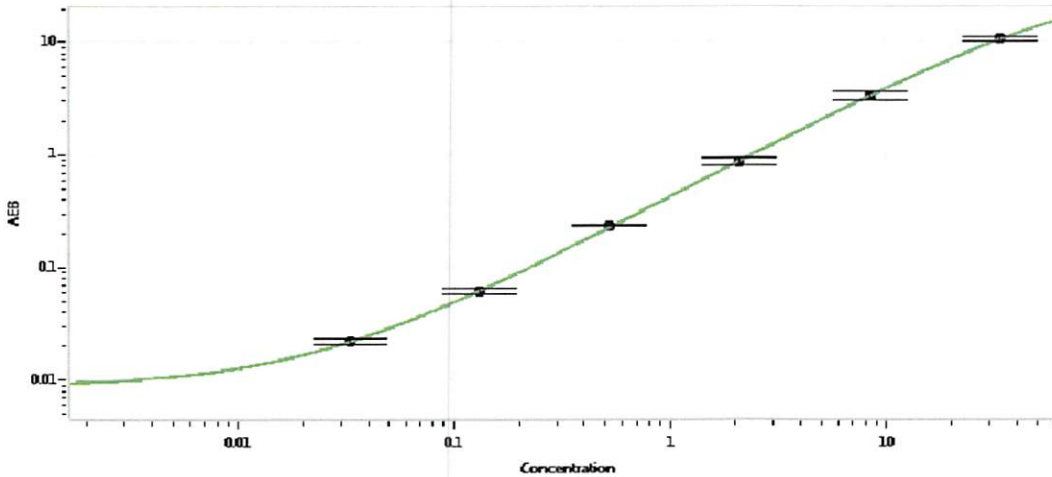
| Component        | Part Number     | Lot Number       |
|------------------|-----------------|------------------|
| Bead Reagent     | 100865          | 819210           |
| Detector Reagent | 100862          | 819111           |
| SBG Reagent      | 100868          | 819021           |
| Sample Diluent   | 100870          | 819022           |
| Calibrators      | 101111          | 81421            |
| Controls         | 100875 & 100876 | 73110            |
| 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          | 73110        | 18.6           | 8.05-24.2          |
| Control 2          | 73110        | 0.343          | 0.137-0.411        |
| Panel 1            | MF-011916-05 | 2.22           | 0.919-2.76         |
| Panel 2            | MF-011916-06 | 16.8           | 5.84-17.5          |

\* 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                         | 0.030 |
| C                         | 0.106 |
| D                         | 0.416 |
| E                         | 1.73  |
| F                         | 6.86  |
| G                         | 26.7  |

Review/Approval

Patrizia Stadler  
Name

Associate Director,  
QC  
Title

*p. Stadler*  
Signature

18 JAN 2019  
Date