

|                 |             |
|-----------------|-------------|
| Product Number: | 101580      |
| Lot Number:     | 501755      |
| Expiration:     | 14-Mar-2020 |
| Platform(s):    | HD-1 & SR-X |

| Component              | Part Number | Lot Number       |
|------------------------|-------------|------------------|
| Bead Reagent           | 101582      | 908803           |
| Detector Reagent       | 101581      | 908510           |
| SBG Reagent            | 101583      | 908407           |
| Sample Diluent         | 103201      | 907309           |
| Calibrator Diluent     | 101585      | 909205           |
| Calibrator Concentrate | 100655      | 909204           |
| RGP Reagent            | 103159      | N/A <sup>1</sup> |

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

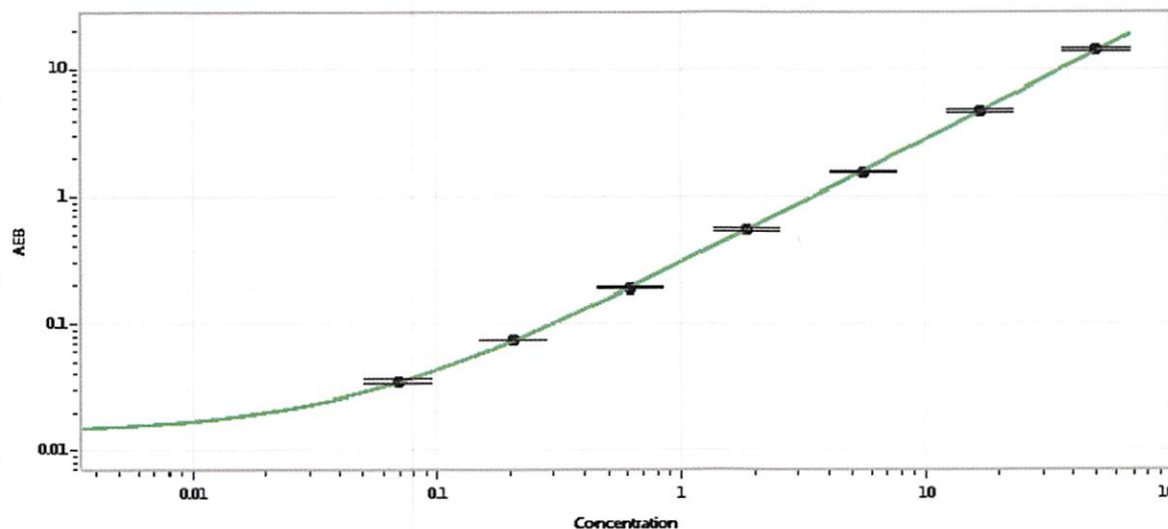


Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Release Materials* |              |                |                    | * Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges. |
|--------------------|--------------|----------------|--------------------|----------------------------------------------------------------------------------------------------------------------|
|                    | Lot          | Result (pg/mL) | Mean Range (pg/mL) |                                                                                                                      |
| Control 1          | MF-042817-01 | 2.90           | 2.05-3.08          |                                                                                                                      |
| Control 2          | MF-051617-02 | 44.9           | 35.8-53.7          |                                                                                                                      |
| Panel 1            | MF-051017-03 | 30.6           | 20.9-32.4          |                                                                                                                      |
| Panel 2            | MF-032217-11 | 2.03           | 1.33-2.15          |                                                                                                                      |
| Panel 3            | MF-033018-06 | 3.21           | 1.68-3.50          |                                                                                                                      |

|                                                 |            |
|-------------------------------------------------|------------|
| Calibrator Concentrate Value Assignment (pg/mL) | 1065 pg/mL |
|-------------------------------------------------|------------|

Calibration Curve



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 0.069 |
| C                         | 0.206 |
| D                         | 0.617 |
| E                         | 1.85  |
| F                         | 5.56  |
| G                         | 16.7  |
| H                         | 50.0  |

|                 |                  |                     |                                 |
|-----------------|------------------|---------------------|---------------------------------|
| Review/Approval | Patrizia Stadler | Assoc. Director, QC | <i>P. Stadler</i> 05 April 2019 |
|                 | Name             | Title               |                                 |