

|                 |             |
|-----------------|-------------|
| Product Number: | 102648      |
| Lot Number:     | 501604      |
| Expiration:     | 24-May-2019 |
| Platform(s):    | HD-1        |

| Component              | Part Number | Lot Number   |
|------------------------|-------------|--------------|
| Bead Stock             | 102641      | 901810       |
| Detector Stock         | 102642      | 901908       |
| Detector Diluent       | 102734      | 901907       |
| Calibrator Concentrate | 102645      | MF-052417-01 |
| Calibrator Diluent     | 102643      | 901014       |
| Sample Diluent         | 102644      | 826227       |

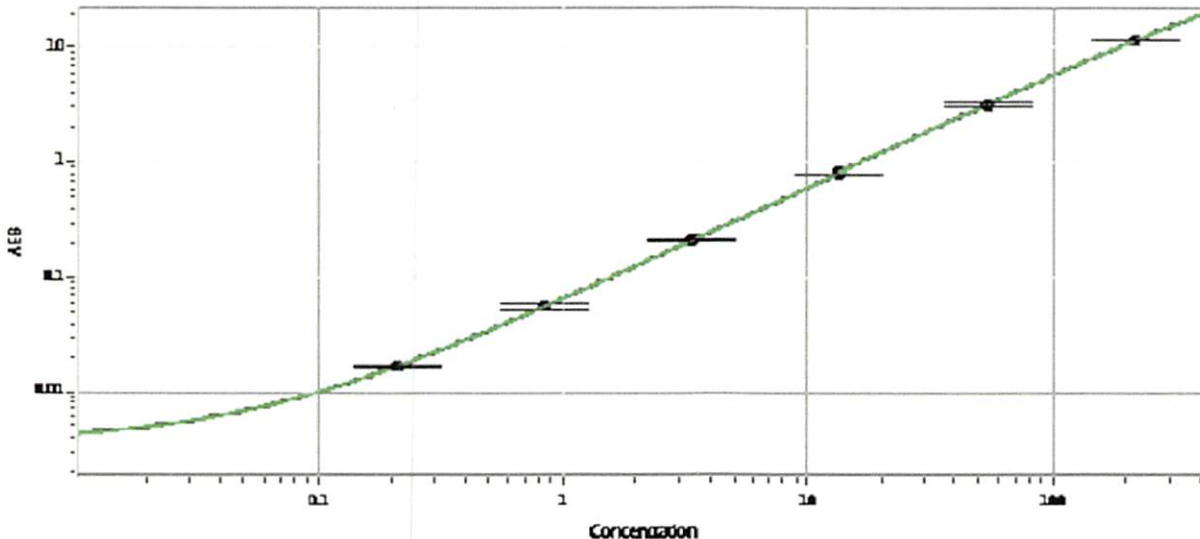


Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Generic Components - Included With Kit But Not Lot Number Specific |        |
|--|--------|
| Discovery Bead Diluent   | 101954 |
| Discovery SBG Diluent  | 101956 |
| Discovery SBG Stock  | 103297 |
| RGP Reagent  | 103159 |

|                                     |             |
|-------------------------------------|-------------|
| Calibrator Concentrate Target Value | 3,000 pg/mL |
|-------------------------------------|-------------|

Calibration Curve



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 0.210 |
| C                         | 0.840 |
| D                         | 3.36  |
| E                         | 13.4  |
| F                         | 53.8  |
| G                         | 215   |

Review/Approval

Patrizia Stadler  
Name

Associate Director, QC  
Title

*P. Stadler* 29 Jan 2019  
Signature/Date