

|                 |             |
|-----------------|-------------|
| Product Number: | 102318      |
| Lot Number:     | 502983      |
| Expiration:     | 17-Mar-2022 |
| Platform(s):    | HD-1/HDX    |



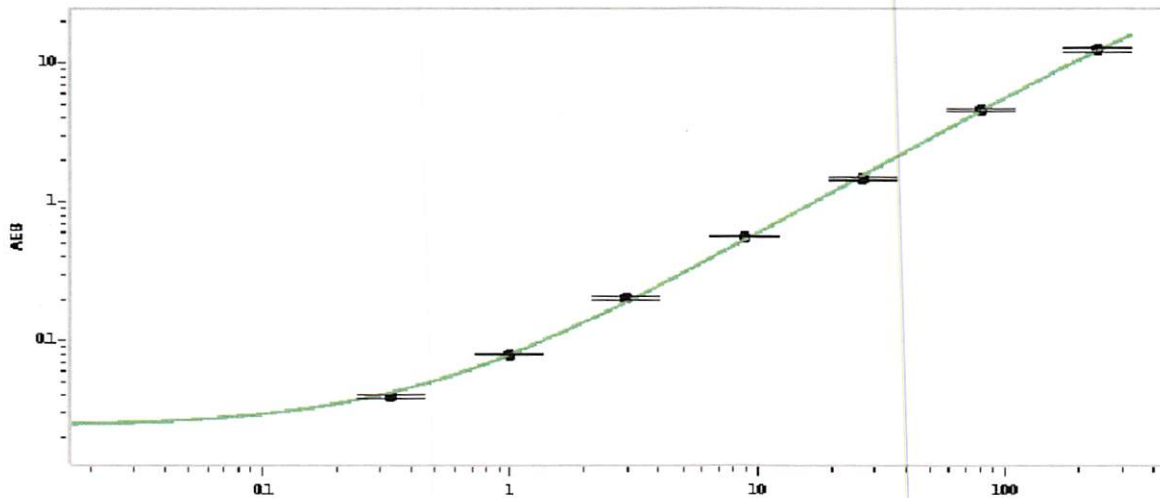
Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Component              | Part Number | Lot Number |
|------------------------|-------------|------------|
| Bead Stock             | 102309      | 107605     |
| Detector Stock         | 102310      | 115808     |
| Calibrator Concentrate | 102314      | 030204     |
| Calibrator Diluent     | 102311      | 107707     |
| Sample Diluent         | 102312      | 107706     |

| Generic Components - Included With Kit But Not Lot Number Specific |        |
|--|--------|
| Discovery Bead Diluent   | 101954 |
| Discovery Detector Diluent   | 101955 |
| Discovery SBG Diluent  | 101956 |
| Discovery SBG Stock  | 103297 |
| RGP Reagent  | 103159 |

|                                     |             |
|-------------------------------------|-------------|
| Calibrator Concentrate Target Value | 4,800 pg/mL |
|-------------------------------------|-------------|

Calibration Curve



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 0.329 |
| C                         | 0.988 |
| D                         | 2.96  |
| E                         | 8.89  |
| F                         | 26.7  |
| G                         | 80.0  |
| H                         | 240   |

Review/Approval

Junpil Park  
Name

Supervisor, QC  
Title

 15 Jun 2021  
Signature/Date