

|                 |             |
|-----------------|-------------|
| Product Number: | 102635      |
| Lot Number:     | 503352      |
| Expiration:     | 10-Nov-2022 |
| Platform(s):    | HD-1/ HD-X  |



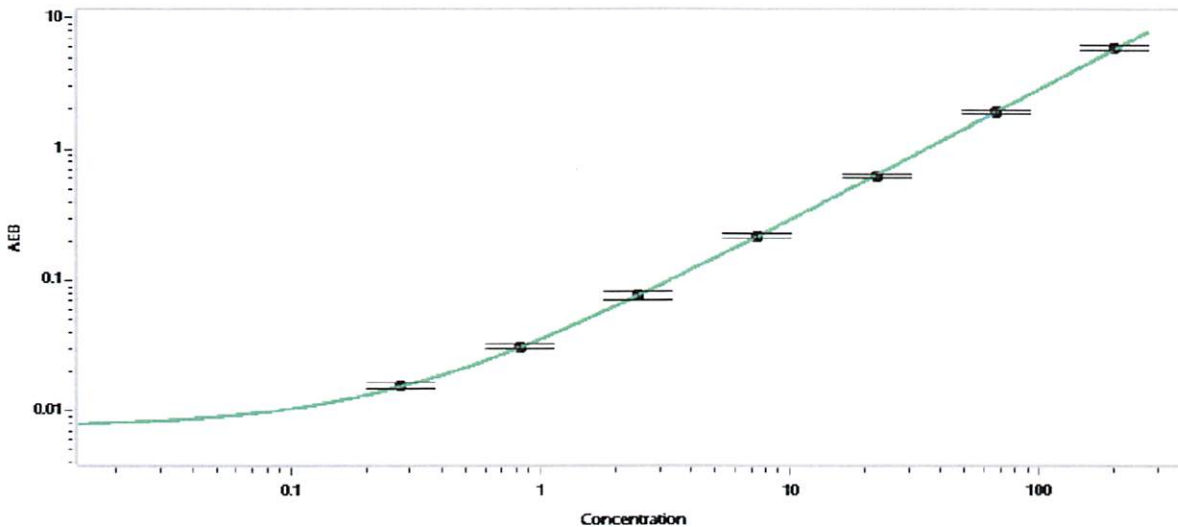
Data below represents results generated on the Simoa® HD-X Analyzer.

| Component              | Part Number | Lot Number |
|------------------------|-------------|------------|
| Bead Stock             | 102628      | 204704     |
| Detector Stock         | 102629      | 215801     |
| Calibrator Concentrate | 102632      | 205503     |
| Calibrator Diluent     | 102630      | 202212     |
| Sample Diluent         | 102631      | 204804     |

| 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 (ng/mL) | 4 |
|---|---|

Calibration Curve



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.00  |
| B                         | 0.274 |
| C                         | 0.823 |
| D                         | 2.47  |
| E                         | 7.41  |
| F                         | 22.2  |
| G                         | 66.7  |
| H                         | 200   |

Review/Approval

Matt Smart Director, QC  
Name Title

M.A. 11/5/2022  
Signature/Date