

| | |
|-----------------|-------------|
| Product Number: | 101984 |
| Lot Number: | 503534 |
| Expiration: | 22-May-2023 |
| Platform(s): | HD-1/HD-X |

| Component | Part Number | Lot Number |
|------------------------|-------------|------------|
| Bead Stock | 101910 | 221001 |
| Detector Stock | 101911 | 221304 |
| Calibrator Concentrate | 101983 | 221407 |
| Calibrator Diluent | 101908 | 221007 |
| Sample Diluent | 101909 | 221006 |

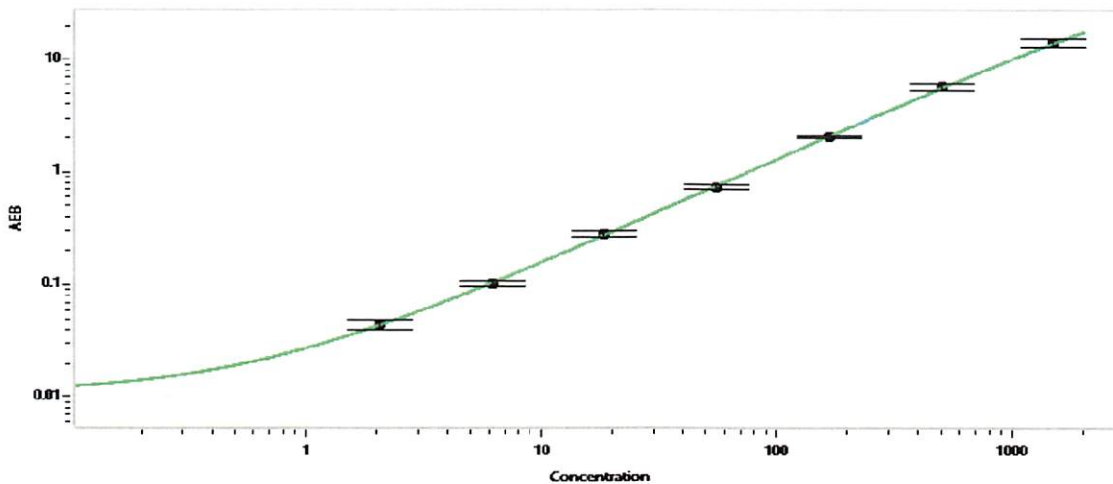


Data below represents results generated on the Simoa® HD-X Analyzer.

| 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) | 30.0 |
|---|------|

Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 2.06 |
| C | 6.17 |
| D | 18.5 |
| E | 55.6 |
| F | 167 |
| G | 500 |
| H | 1500 |

Review/Approval: Sonia Khillan Name, QC Manager Title, Sonia Khillan 29 AUG 2022 Signature/Date