

| | |
|------------------------|--------------------|
| Product Number: | 102336 |
| Lot Number: | 504233 |
| Expiration: | 7-Nov-2025 |
| Platform(s): | SR-X and HD-1/HD-X |



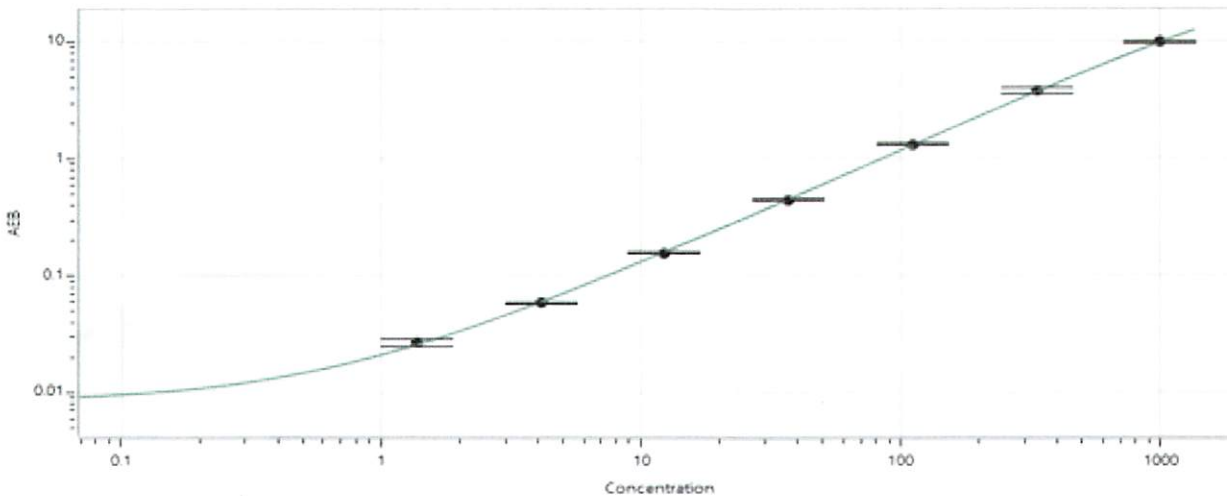
| Component | Part Number | Lot Number |
|------------------------|-------------|------------|
| Bead Stock | 102330 | 412801 |
| Detector Stock | 102331 | 416906 |
| Calibrator Concentrate | 102335 | 414303 |
| Calibrator Diluent | 102332 | 412842 |
| Sample Diluent | 102333 | 412803 |

| 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 |

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

| | |
|--|----|
| Calibrator Concentrate Target Value (ng/mL) | 20 |
|--|----|

Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 1.37 |
| C | 4.12 |
| D | 12.3 |
| E | 37.0 |
| F | 111 |
| G | 333 |
| H | 1000 |

Review/Approval

Cody Carlson
Name

QC Manager
Title

[Signature] 21 JUN 2024
Signature/Date