

| | |
|------------------------|----------|
| Product Number: | 102233 |
| Lot Number: | 501076 |
| Expiration: | 5-Jun-18 |
| Platform(s): | HD-1 |

| Component | Part Number | Lot Number |
|------------------------|-------------|--------------|
| Bead Stock | 102223 | MF-010518-14 |
| Detector Stock | 102224 | MF-120517-06 |
| Calibrator Concentrate | 102229 | MF-112116-01 |
| Calibrator Diluent | 102225 | MF-120517-07 |
| Sample Diluent | 102226 | MF-120517-08 |

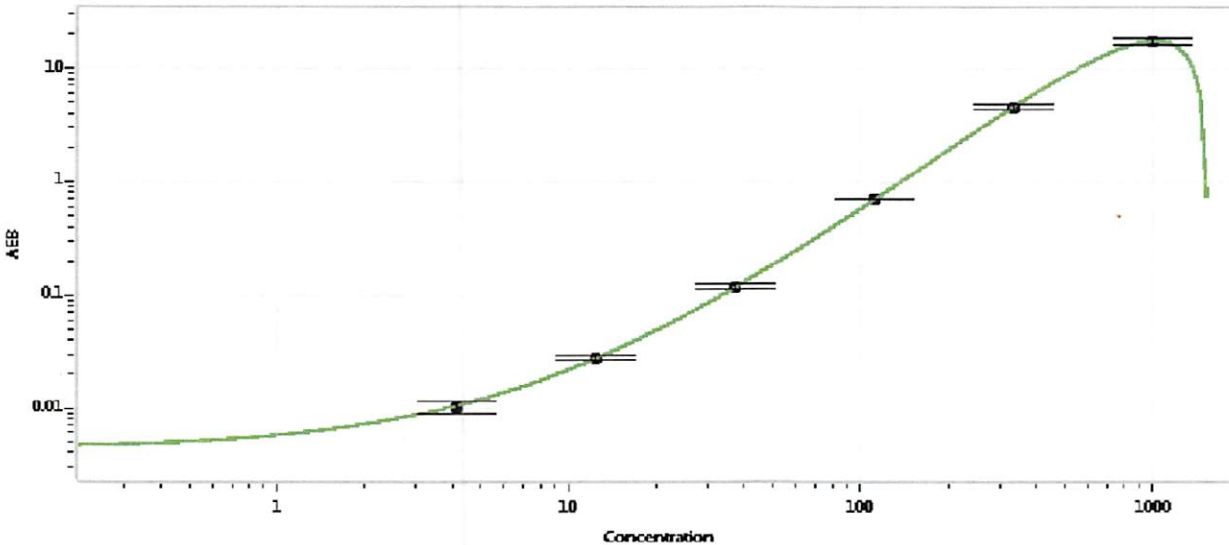


Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Generic Components - Included With Kit But Not Lot Number Specific | |
|--|--------|
| Discovery Bead Diluent | 101954 |
| Discovery Detector Diluent | 101955 |
| Discovery SBG Diluent | 101956 |
| Discovery SBG Concentrate | 101957 |
| RGP Reagent | 103159 |

| | |
|--|----------|
| Alpha-Synuclein Calibrator Concentrate Target Value | 20 ng/mL |
|--|----------|

Calibration Curve



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

Review/Approval: Sivasankari Murugan Name, QC Supervisor Title, [Signature] Signature/Date, 22 Jan 2018 Date