

| | |
|-----------------|------------|
| Product Number: | 102073 |
| Lot Number: | 502485 |
| Expiration: | 2-May-2021 |
| Platform(s): | HD-1/HD-X |

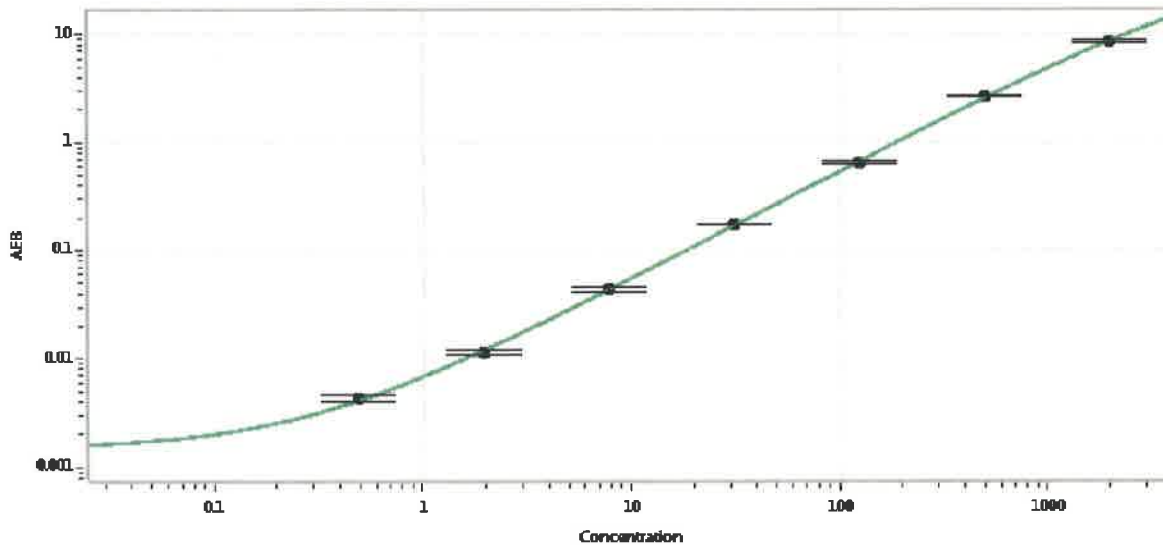


| Component | Part Number | Lot Number |
|------------------------|-------------|------------|
| Bead Stock | 102065 | 018906 |
| Detector Stock | 102066 | 021103 |
| Calibrator Concentrate | 102069 | 019101 |
| Calibrator Diluent | 102067 | 019006 |
| Sample Diluent | 102068 | 019005 |

| 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-1 Analyzer.

| | |
|-------------------------------------|----------|
| Calibrator Concentrate Target Value | 40 ng/mL |
|-------------------------------------|----------|



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.488 |
| C | 1.95 |
| D | 7.81 |
| E | 31.3 |
| F | 125 |
| G | 500 |
| H | 2000 |

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 06 Aug 2020

Name

Title

Signature/Date