

| | |
|-----------------|--------------------|
| Product Number: | 103714 |
| Lot Number: | 503199 |
| Expiration: | 29-Oct-2022 |
| Platform(s): | HD-1 / HD-X & SR-X |



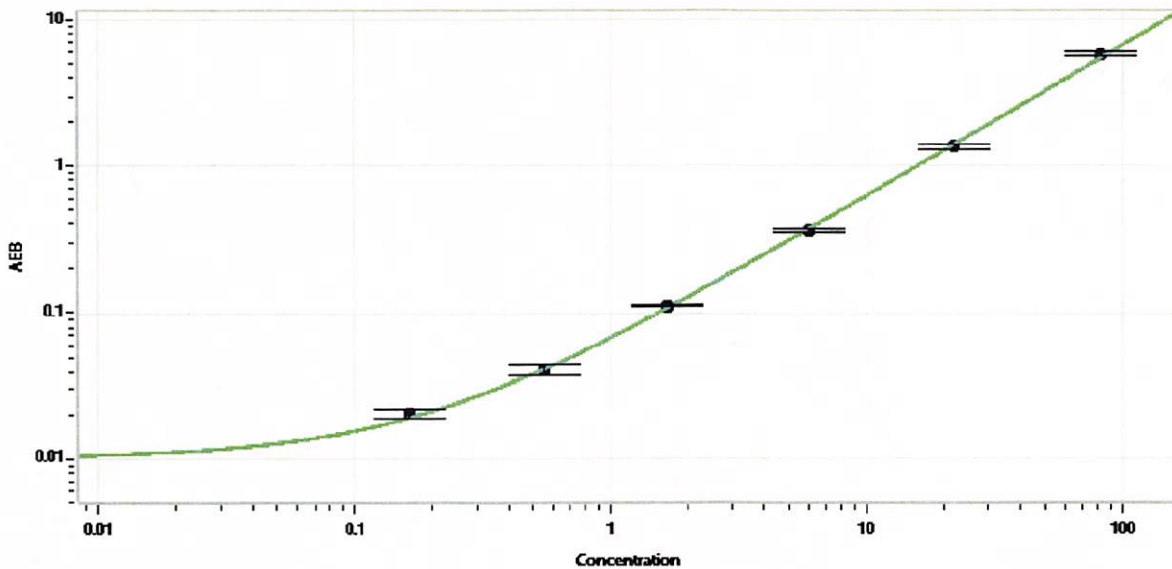
Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Component | Part Number | Lot Number |
|----------------------|-------------|------------------|
| Bead Reagent | 103698 | 130207 |
| Detector Reagent | 103700 | 130513 |
| SBG Reagent | 103702 | 130403 |
| Sample Diluent | 103704 | 130405 |
| Reference Calibrator | 103711 | 134201 |
| Control 1 | 103712 | 134202 |
| Control 2 | 103713 | 202510 |
| RGP Reagent | 103159 | N/A ¹ |

¹ RGP is not Kit Lot Specific

| Release Materials* | | | | * Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges. |
|--------------------|--------|----------------|--------------------|--|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) | |
| Control 1 | 134202 | 3.82 | 3.03-4.91 | |
| Control 2 | 202510 | 105 | 81.7-123 | |
| Panel 1 | 013903 | 0.668 | 0.258-1.11 | |
| Panel 2 | 013903 | 1.98 | 0.841-2.44 | |
| Panel 3 | 013903 | 3.71 | 1.79-4.31 | |

Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.165 |
| C | 0.552 |
| D | 1.67 |
| E | 5.99 |
| F | 22.0 |
| G | 82.0 |

Review/Approval: Shivani Goel Name QC Title Shivani Goel Signature/Date 11 Feb 2022