

| | |
|-----------------|-------------|
| Product Number: | 103186 |
| Lot Number: | 503168 |
| Expiration: | 27-Aug-2022 |
| Platform(s): | HD-1/ HD-X |



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

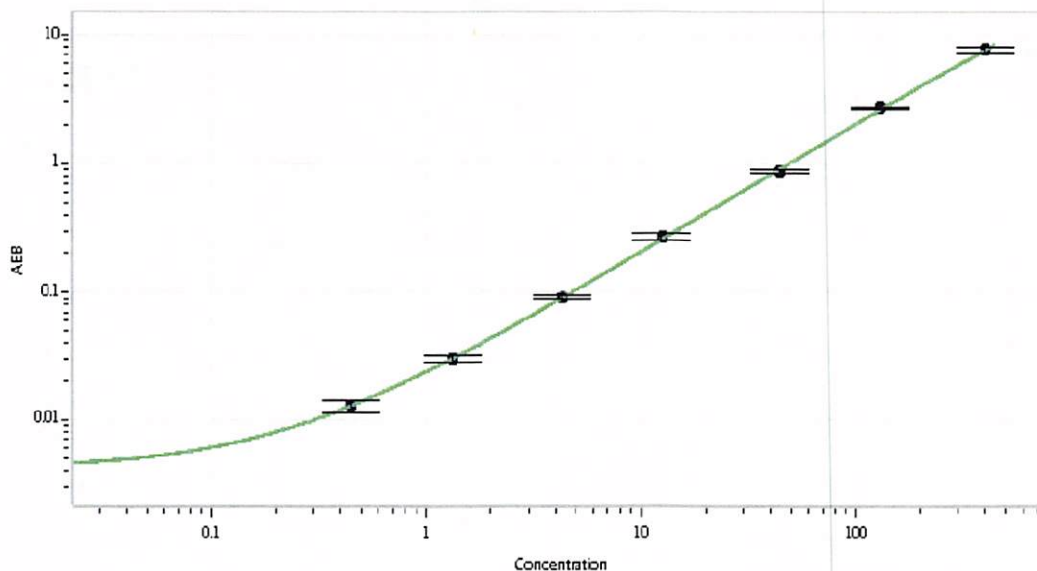
| Component | Part Number | Lot Number |
|------------------|-------------|------------------|
| Bead Reagent | 102246 | 124505 |
| Detector Reagent | 102248 | 123907 |
| SBG Reagent | 102250 | 127405 |
| Sample Diluent | 102252 | 123909 |
| Calibrators | 102255 | 133605 |
| Control 1 | 102256 | 133606 |
| Control 2 | 102257 | 133606 |
| RGP Reagent | 103159 | N/A ¹ |

¹ RGP is not Kit Lot Specific

| Release Materials* | | | |
|--------------------|--------|----------------|--------------------|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 133606 | 4.32 | 3.12 - 5.20 |
| Control 2 | 133606 | 264 | 202 - 303 |
| Panel 1 | 008501 | 13.7 | 8.59 - 20.5 |
| Panel 2 | 008501 | 10.7 | 7.39 - 15.1 |
| Panel 3 | 008501 | 63.4 | 43.8 - 91.8 |
| Panel 4 | 009401 | 193 | 128 - 302 |

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 0.447 |
| C | 1.33 |
| D | 4.33 |
| E | 12.5 |
| F | 44.2 |
| G | 129 |
| H | 400 |

Review/Approval Shivani Goel QC Shivani Goel 30 Dec 2021
Name Title Signature/Date