

| | |
|-----------------|-------------|
| Product Number: | 103186 |
| Lot Number: | 503215 |
| Expiration: | 27-Jun-2022 |
| Platform(s): | HD-1/ HD-X |



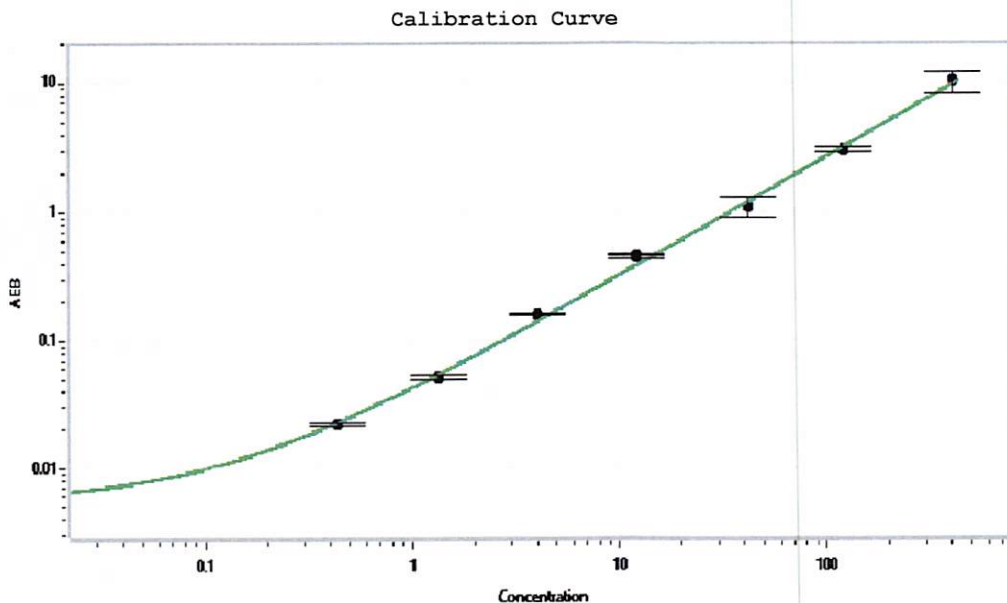
Data below represents results generated on the Simoa HD-1® Analyzer.

| Component | Part Number | Lot Number |
|------------------|-------------|------------------|
| Bead Reagent | 102246 | 117801 |
| Detector Reagent | 102248 | 117803 |
| SBG Reagent | 102250 | 117808 |
| Sample Diluent | 102252 | 117914 |
| Calibrators | 102255 | 130504 |
| Control 1 | 102256 | 130505 |
| Control 2 | 102257 | 130505 |
| 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 | 130505 | 3.62 | 2.99 - 4.57 |
| Control 2 | 130505 | 203 | 111 - 225 |
| Panel 1 | 008501 | 13.8 | 8.59 - 20.5 |
| Panel 2 | 008501 | 10.8 | 7.39 - 15.1 |
| Panel 3 | 008501 | 71.2 | 43.8 - 91.8 |
| Panel 4 | 009401 | 275 | 128.4 - 302.1 |

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



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.439 |
| C | 1.35 |
| D | 4.02 |
| E | 12.0 |
| F | 41.6 |
| G | 120 |
| H | 407 |

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature]
Signature/Date

18 Nov 2021