

| | |
|-----------------|------------------|
| Product Number: | 102732 |
| Lot Number: | 503123 |
| Expiration: | 2-Aug-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 | 102722 | 124210 |
| Detector Reagent | 102724 | 124303 |
| SBG Reagent | 102726 | 124205 |
| Sample Diluent | 102728 | 124008 |
| Calibrator Diluent | 102731 | 124007 |
| Calibrator Concentrate | 100658 | 124304 |
| RGP Reagent | 103159 | N/A ¹ |

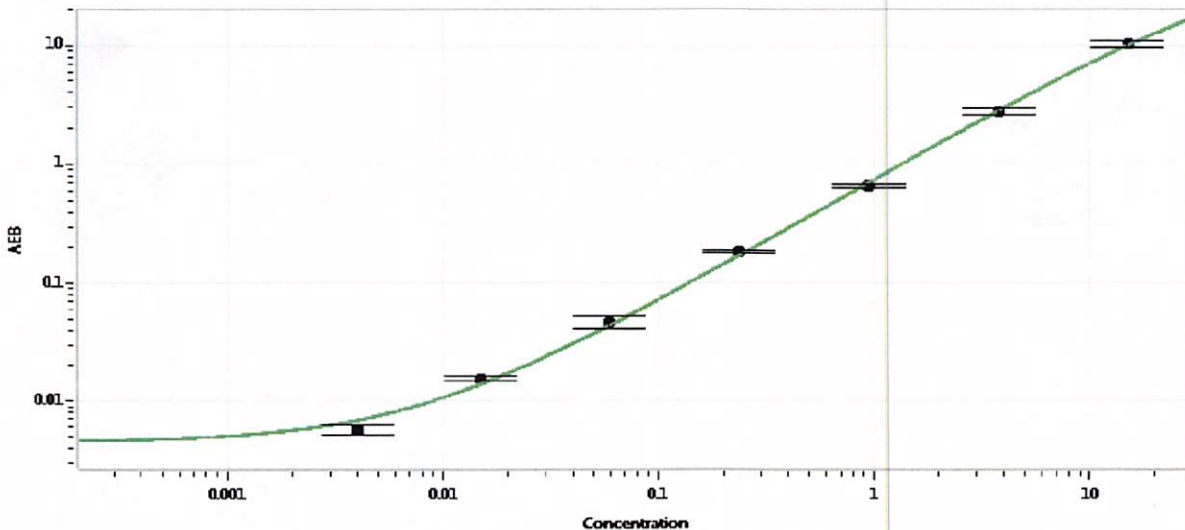
¹ RGP is not Kit Lot Specific

| Release Materials* | | | |
|--------------------|--------|----------------|--------------------|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 114108 | 0.110 | 0.067 - 0.159 |
| Control 2 | 114108 | 2.40 | 1.91 - 3.81 |
| Panel 2 | 112714 | 0.034 | 0.010 - 0.049 |
| Panel 3 | 112714 | 1.46 | 1.02 - 1.87 |
| Panel 4 | 112714 | 5.49 | 2.78 - 7.70 |

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

| | |
|---|-----|
| Calibrator Concentrate Value Assignment (pg/mL) | 422 |
|---|-----|

Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.004 |
| C | 0.015 |
| D | 0.059 |
| E | 0.234 |
| F | 0.938 |
| G | 3.75 |
| H | 15.0 |

Review/Approval

Junpil Park

QC Supervisor

[Signature] 09sep2021

Name

Title

Signature/Date