

| | |
|-----------------|------------------|
| Product Number: | 102732 |
| Lot Number: | 502378 |
| Expiration: | 16-Apr-2021 |
| 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 | 012807 |
| Detector Reagent | 102724 | 012804 |
| SBG Reagent | 102726 | 012805 |
| Sample Diluent | 102728 | 012806 |
| Calibrator Diluent | 102731 | 010715 |
| Calibrator Concentrate | 100658 | 013605 |
| RGP Reagent | 103159 | N/A ¹ |

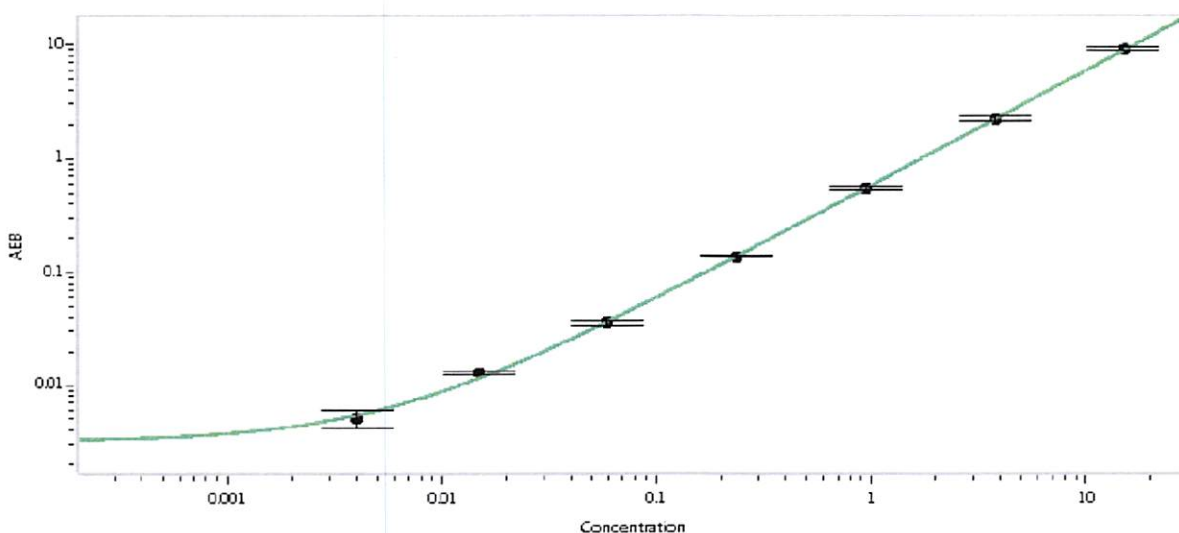
¹ RGP is not Kit Lot Specific

| Release Materials* | | | |
|--------------------|--------|----------------|--------------------|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 909502 | 0.124 | 0.095-0.162 |
| Control 2 | 909502 | 3.25 | 2.88-4.52 |
| Panel 1 | 911401 | 0.026 | 0.013-0.050 |
| Panel 3 | 911403 | 1.53 | 1.22-2.02 |
| Panel 4 | 911404 | 5.15 | 3.95-7.43 |

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

| | |
|---|-----|
| Calibrator Concentrate Value Assignment (pg/mL) | 497 |
|---|-----|

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: Shivani Goel QC Manager Shivani Goel 28 May 2020
Name Title Signature/Date