

| | |
|-----------------|-------------|
| Product Number: | 102732 |
| Lot Number: | 501204 |
| Expiration: | 26-Aug-2018 |
| Platform(s): | HD-1 & SR-X |



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

| Component | Part Number | Lot Number |
|------------------------|-------------|------------------|
| Bead Reagent | 102722 | 80953 |
| Detector Reagent | 102724 | 80932 |
| SBG Reagent | 102726 | 80933 |
| Sample Diluent | 102728 | 80573 |
| Calibrator Diluent | 102731 | 80943 |
| Calibrator Concentrate | 100658 | 73217 |
| RGP Reagent | 103159 | N/A ¹ |

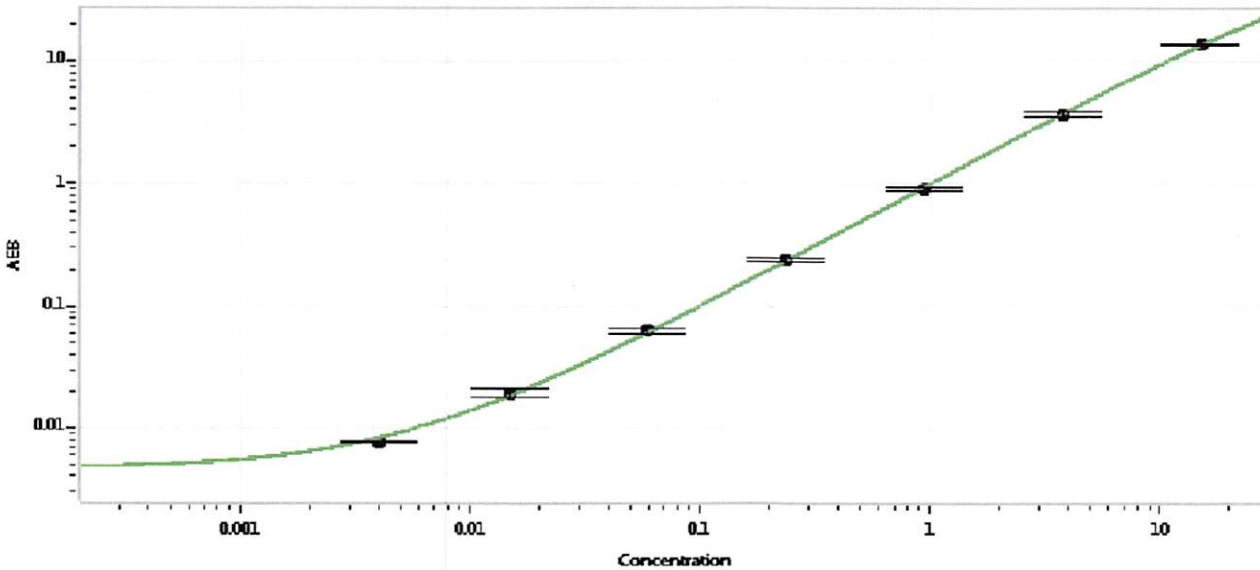
¹ RGP is not Kit Lot Specific

| Release Materials* | | | |
|--------------------|--------------|----------------|--------------------|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | MF-061217-01 | 0.201 | 0.134-0.289 |
| Control 2 | MF-061217-01 | 3.72 | 2.44-6.35 |
| Panel 1 | MF-061317-03 | 0.0735 | 0.0371-0.0876 |
| Panel 2 | MF-061317-04 | 0.335 | 0.191-0.490 |
| Panel 3 | MF-061317-05 | 4.84 | 3.18-7.89 |

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

| | |
|---|-----|
| Calibrator Concentrate Value Assignment (pg/mL) | 194 |
|---|-----|

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

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date 12 MAR 2018