

| | |
|-----------------|------------|
| Product Number: | 100196 |
| Lot Number: | 501339 |
| Expiration: | 7-Nov-2018 |
| Platform(s): | HD-1 |



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

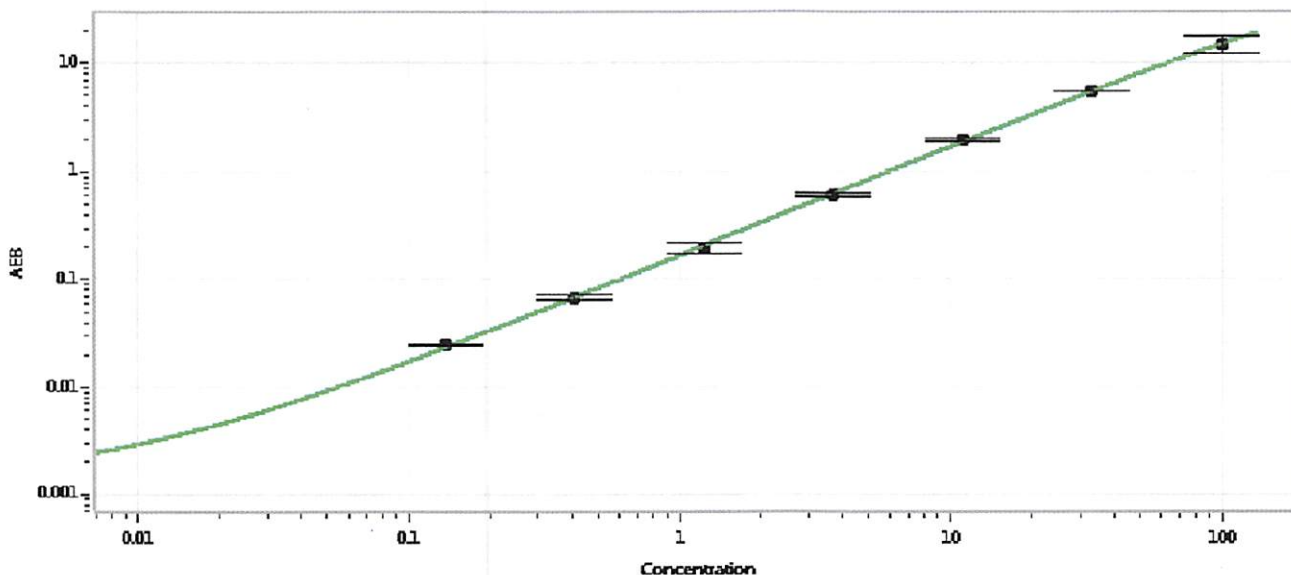
| Component | Part Number | Lot Number |
|------------------------|-------------|------------------|
| Bead Reagent | 100391 | 818410 |
| Detector Reagent | 100315 | 818708 |
| SBG Reagent | 100316 | 818000 |
| Sample Diluent | 100317 | 819709 |
| Calibrator Diluent | 100665 | 81273 |
| Calibrator Concentrate | 100664 | 72701 |
| RGP Reagent | 103159 | N/A ¹ |

¹ RGP is not Kit Lot Specific

| Release Materials* | | | | * Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges. |
|--------------------|--------------|----------------|--------------------|--|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) | |
| Control 1 | 52521 | 98.5 | 75.6-118 | |
| Control 2 | 52521 | 10.1 | 7.72-11.6 | |
| Panel 1 | MF-012116-05 | 1.34 | 0.871-1.73 | |
| Panel 2 | MF-090815-18 | 7.39 | 4.45-9.13 | |
| Panel 3 | MF-090815-19 | 72.3 | 53.9-94.9 | |

| | |
|---|------|
| Calibrator Concentrate Value Assignment (pg/mL) | 2154 |
|---|------|

Calibration Curve



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 0.137 |
| C | 0.412 |
| D | 1.23 |
| E | 3.70 |
| F | 11.1 |
| G | 33.3 |
| H | 100 |

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 23/11/2018
Signature/Date