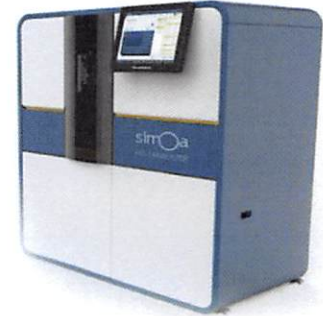


|                 |             |
|-----------------|-------------|
| Product Number: | 101160      |
| Lot Number:     | 501437      |
| Expiration:     | 12-Dec-2018 |
| Platform(s):    | HD-1 & SR-X |



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

| Component            | Part Number     | Lot Number       |
|----------------------|-----------------|------------------|
| Bead Reagent †       | 101263          | 81631            |
| Detector Reagent     | 101264          | 820604           |
| SBG Reagent          | 101296          | 823212           |
| Sample Diluent       | 103314          | 818008           |
| Calibrators          | 101299          | 81720   826802   |
| Controls             | 101302 & 101316 | 81721            |
| SBG Diluent (SR-X)†† | 103158          | N/A <sup>1</sup> |
| RGP Reagent          | 103159          | N/A <sup>1</sup> |

† Follow the Simoa™ Cytokine 3-Plex A Advantage Kit Instructions (KI-0011 - HD1/KI-0101 - SR-X) to prepare the Bead Reagent.  
†† Specific for use on the SR-X platform - refer to the Simoa™ Cytokine 3-Plex A SR-X Kit Instructions (KI-0101).

<sup>1</sup> Reagents are not Kit Lot Specific

**Release Materials**

**IL-6**

|           | Lot          | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 81721        | 0.929          | 0.786-1.18         |
| Control 2 | 81721        | 26.4           | 21.6-32.3          |
| Panel 2   | MF-062818-02 | 0.919          | 0.765-1.15         |
| Panel 3   | MF-062818-03 | 1.80           | 0.976-2.21         |
| Panel 4   | MF-062818-04 | 13.2           | 9.75-15.2          |

**IL-10**

|           | Lot          | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 81721        | 0.702          | 0.623-0.935        |
| Control 2 | 81721        | 13.4           | 10.9-22.8          |
| Panel 2   | MF-062818-02 | 11.4           | 9.90-18.6          |
| Panel 3   | MF-062818-03 | 0.588          | 0.479-0.935        |
| Panel 4   | MF-062818-04 | 1.23           | 1.10-1.64          |

**TNFα**

|           | Lot          | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 81721        | 1.99           | 1.65-2.48          |
| Control 2 | 81721        | 49.9           | 42.8-64.2          |
| Panel 2   | MF-062818-02 | 1.89           | 1.19-3.18          |
| Panel 3   | MF-062818-03 | 39.5           | 29.9-61.0          |
| Panel 4   | MF-062818-04 | 6.98           | 5.98-9.43          |

Ranges shown are generated internally for new lot release only.

Customer should generate their own control ranges.

Review/Approval

Sivasankari Murugan

QC Supervisor

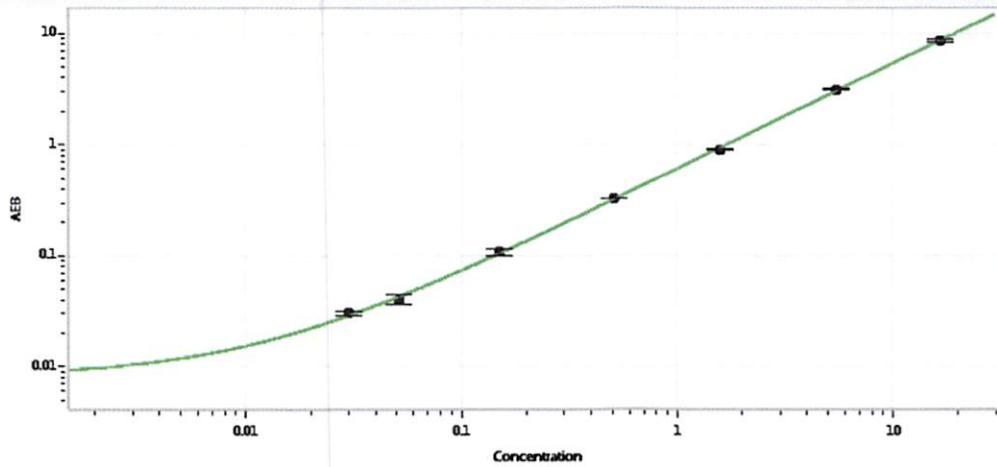
*[Signature]* 050072018

Name

Title

Signature/Date

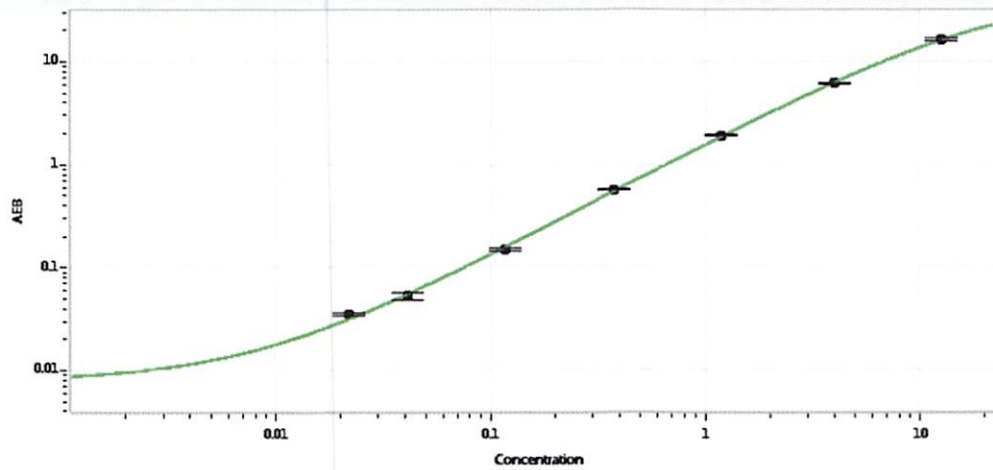
IL-6 Calibration Curve



Calibrator Levels  
(pg/mL)

|   |       |
|---|-------|
| A | 0.000 |
| B | 0.030 |
| C | 0.051 |
| D | 0.150 |
| E | 0.511 |
| F | 1.58  |
| G | 5.50  |
| H | 16.6  |

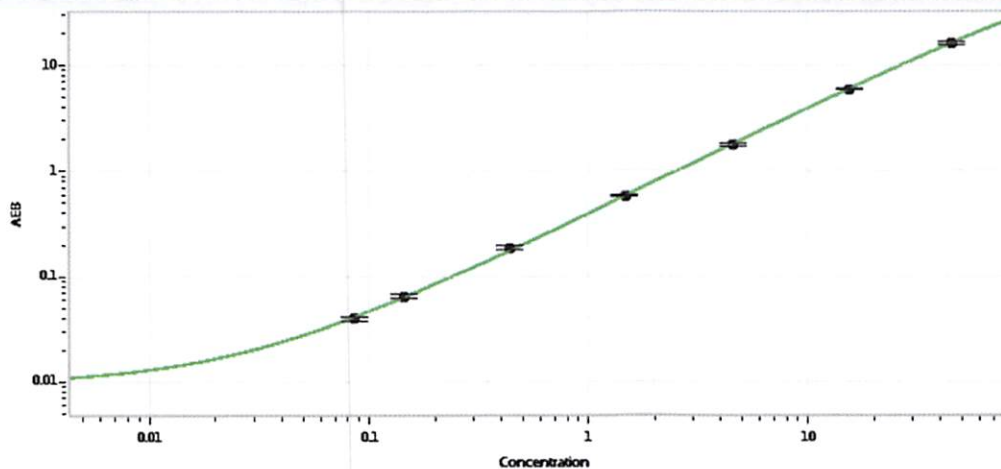
IL-10 Calibration Curve



Calibrator Levels  
(pg/mL)

|   |       |
|---|-------|
| A | 0.000 |
| B | 0.022 |
| C | 0.041 |
| D | 0.117 |
| E | 0.375 |
| F | 1.20  |
| G | 4.02  |
| H | 12.6  |

TNFα Calibration Curve



Calibrator Levels  
(pg/mL)

|   |       |
|---|-------|
| A | 0.000 |
| B | 0.086 |
| C | 0.145 |
| D | 0.440 |
| E | 1.47  |
| F | 4.56  |
| G | 15.4  |
| H | 44.9  |