

| | |
|------------------------|-----------------|
| Product Number: | 101160 |
| Lot Number: | 501951 |
| Expiration: | 13-Jun-2020 |
| 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 † | 101263 | 923409 |
| Detector Reagent | 101264 | 923405 |
| SBG Reagent | 101296 | 920307 |
| Sample Diluent | 103329 | 919810 |
| Calibrators | 103228 | 921801 |
| Controls | 103229-103230 | 921802 |
| SBG Diluent (SR-X)†† | 103158 | N/A ¹ |
| RGP Reagent | 103159 | N/A ¹ |

† 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).

¹ Reagents are not Kit Lot Specific

Release Materials

IL-6

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 921802 | 0.735 | 0.600-0.900 |
| Control 2 | 921802 | 20.8 | 16.8-25.2 |
| Panel 2 | MF-062818-02 | 0.992 | 0.777-1.17 |
| Panel 3 | MF-062818-03 | 1.62 | 1.37-2.06 |
| Panel 4 | MF-062818-04 | 14.4 | 10.4-15.7 |

IL-10

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 921802 | 0.466 | 0.379-0.568 |
| Control 2 | 921802 | 10.6 | 6.85-12.5 |
| Panel 2 | MF-062818-02 | 15.1 | 10.0-15.8 |
| Panel 3 | MF-062818-03 | 1.08 | 0.302-1.18 |
| Panel 4 | MF-062818-04 | 1.65 | 0.897-1.70 |

TNFα

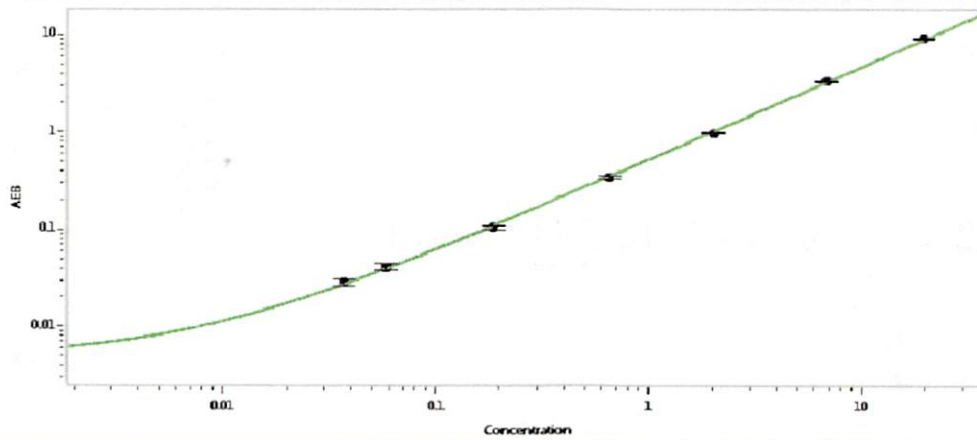
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 921802 | 2.59 | 2.03-3.04 |
| Control 2 | 921802 | 63.5 | 51.0-76.5 |
| Panel 2 | MF-062818-02 | 1.98 | 1.54-2.41 |
| Panel 3 | MF-062818-03 | 39.8 | 31.9-47.8 |
| Panel 4 | MF-062818-04 | 7.17 | 5.51-8.26 |

Ranges shown are generated internally for new lot release only.

Customer should generate their own control ranges.

| | | | | |
|------------------------|------------------|---------------------|-------------------|--------------|
| Review/Approval | Patrizia Stadler | Assoc. Director, QC | <i>P. Stadler</i> | 05 Sept 2019 |
| | Name | Title | Signature | Date |

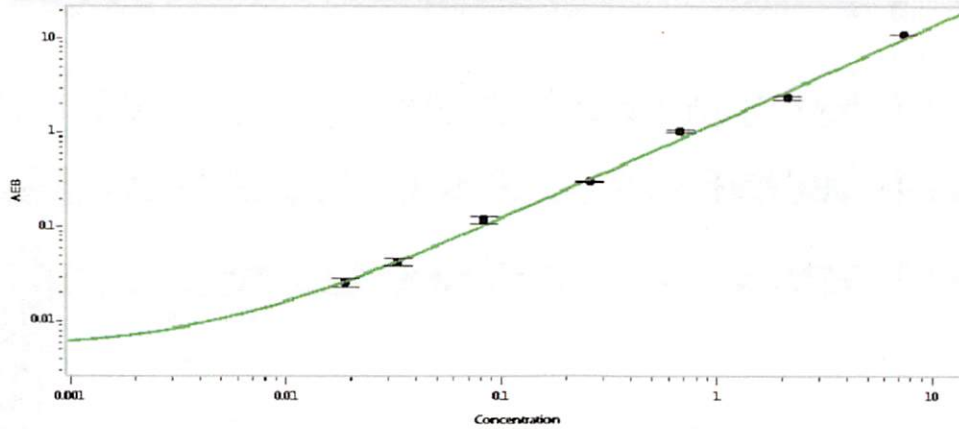
IL-6 Calibration Curve



**Calibrator Levels
(pg/mL)**

| | |
|---|-------|
| A | 0.000 |
| B | 0.037 |
| C | 0.059 |
| D | 0.186 |
| E | 0.653 |
| F | 2.01 |
| G | 6.80 |
| H | 19.6 |

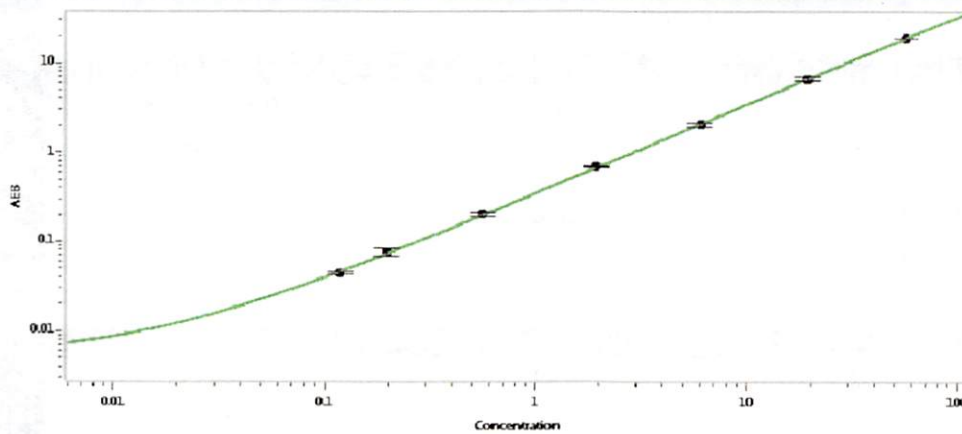
IL-10 Calibration Curve



**Calibrator Levels
(pg/mL)**

| | |
|---|-------|
| A | 0.000 |
| B | 0.019 |
| C | 0.033 |
| D | 0.083 |
| E | 0.258 |
| F | 0.677 |
| G | 2.12 |
| H | 7.33 |

TNFα Calibration Curve



**Calibrator Levels
(pg/mL)**

| | |
|---|-------|
| A | 0.000 |
| B | 0.118 |
| C | 0.197 |
| D | 0.563 |
| E | 1.93 |
| F | 6.04 |
| G | 19.6 |
| H | 56.5 |