

| | |
|-----------------|-------------|
| Product Number: | 101319 |
| Lot Number: | 502000 |
| Expiration: | 17-Jul-2020 |
| Platform(s): | HD-1/HD-x |



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

| Component | Part Number | Lot Number |
|------------------|---------------|------------------|
| Bead Reagent † | 101265 | 922401 |
| Detector Reagent | 101266 | 923306 |
| SBG Reagent | 101297 | 920308 |
| Sample Diluent | 103330 | 919811 |
| Calibrators | 103228 | 921801 |
| Controls | 103229-103230 | 921802 |
| RGP Reagent | 103159 | N/A ¹ |

† Follow the Simoa™ Cytokine 3-Plex B Advantage Kit Instructions (KI-0012) to prepare the Bead Reagent.

¹ Reagents are not Kit Lot Specific

Release Materials

IL-6

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 921802 | 0.711 | 0.579-0.868 |
| Control 2 | 921802 | 20.5 | 16.9-25.3 |
| Panel 2 | MF-062818-02 | 0.986 | 0.763-1.14 |
| Panel 3 | MF-062818-03 | 1.67 | 1.28-1.91 |
| Panel 4 | MF-062818-04 | 14.7 | 11.0-16.5 |

IL-17A

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 921802 | 0.523 | 0.426-0.639 |
| Control 2 | 921802 | 11.0 | 9.39-14.1 |
| Panel 2 | MF-062818-02 | 18.6 | 13.9-20.8 |
| Panel 3 | MF-062818-03 | 1.51 | 1.18-1.76 |
| Panel 4 | MF-062818-04 | 1.95 | 1.59-2.38 |

TNFα

| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
|-----------|--------------|----------------|--------------------|
| Control 1 | 921802 | 2.28 | 1.84-2.76 |
| Control 2 | 921802 | 59.3 | 50.0-75.0 |
| Panel 2 | MF-062818-02 | 1.64 | 1.40-2.55 |
| Panel 3 | MF-062818-03 | 41.1 | 32.6-51.5 |
| Panel 4 | MF-062818-04 | 6.39 | 5.49-8.23 |

Ranges shown are generated internally for new lot release only.

Customer should generate their own control ranges.

Review/Approval

Daniel Ritchie Senior QC Analyst

Name

Title

[Signature] / 22 Aug 2019
Signature/Date

*Signing on behalf of Patrizia Stadler