



# CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07039

**Product Description:** Cytokine Panel 1 (human) Control Pack

**Catalog Number:** C4050-1

**Lot Number:** CP000039

**Expiration Date:** 30 APR 2020

## Concentration of Materials Tested:

| Units = pg/mL                      | Assay     |                  |         |         |                 |          |          |           |                 |         |
|------------------------------------|-----------|------------------|---------|---------|-----------------|----------|----------|-----------|-----------------|---------|
|                                    | hu GM-CSF | hu IL-1 $\alpha$ | hu IL-5 | hu IL-7 | hu IL-12/23 p40 | hu IL-15 | hu IL-16 | hu IL-17A | hu TNF- $\beta$ | hu VEGF |
| Reconstituted Control 1 (A00C0325) | 362       | 167              | 371     | 241     | 1290            | 241      | 1160     | 2590      | 196             | 373     |
| Reconstituted Control 2 (A00C0326) | 55.8      | 57.8             | 123     | 42.0    | 208             | 36.5     | 328      | 327       | 31.7            | 127     |
| Reconstituted Control 3 (A00C0393) | 9.65      | 19.3             | 32.8    | 7.49    | 35.1            | 8.01     | 110      | 57.1      | 5.71            | 49.0    |

**Reconstitution Volume:** 250  $\mu$ L

Control concentrations were assigned across three kit lots.

| Cytokine Panel 1 (human) Control Pack Components |            |                     |                 |
|--|------------|---------------------|-----------------|
| Description                                      | Lot Number | Storage Temperature | Expiration Date |
| Cytokine Panel 1 (human) Control 1               | A00C0325   | 2-8°C               | 30 APR 2020     |
| Cytokine Panel 1 (human) Control 2               | A00C0326   | 2-8°C               | 31 MAY 2020     |
| Cytokine Panel 1 (human) Control 3               | A00C0393   | 2-8°C               | 30 APR 2021     |

**Statement:** The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

|                        | Name                | Function          | Signature                  | Date        |
|------------------------|---------------------|-------------------|----------------------------|-------------|
| <b>Review/Approval</b> | Vanessa Castleberry | Quality Assurance | <i>Vanessa Castleberry</i> | 21 JUN 2016 |