



### CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-04462

**Product Description:** Proinflammatory Panel 1 (human) Control Pack  
**Catalog Number:** C4049-1  
**Lot Number:** CP000016  
**Expiration Date:** 31 MAY 2019

**Concentration of Materials Tested:**

| Units = pg/mL                      | Assay            |                 |         |         |         |         |          |             |          |                  |
|------------------------------------|------------------|-----------------|---------|---------|---------|---------|----------|-------------|----------|------------------|
|                                    | hu IFN- $\gamma$ | hu IL-1 $\beta$ | hu IL-2 | hu IL-4 | hu IL-6 | hu IL-8 | hu IL-10 | hu IL-12p70 | hu IL-13 | hu TNF- $\alpha$ |
| Reconstituted Control 1 (A00C0247) | 401              | 136             | 287     | 55.6    | 246     | 157     | 98.3     | 109         | 152      | 60.5             |
| Reconstituted Control 2 (A00C0248) | 106              | 32.8            | 70.5    | 13.2    | 56.8    | 37.6    | 24.3     | 23.6        | 33.2     | 12.4             |
| Reconstituted Control 3 (A00C0249) | 39.9             | 7.71            | 17.0    | 3.19    | 11.8    | 8.52    | 5.44     | 7.26        | 15.2     | 3.88             |

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

Control concentrations were assigned across three kit lots.

| Proinflammatory Panel 1 (human) Control Pack Components |            |                     |                 |
|---|------------|---------------------|-----------------|
| Description   | Lot Number | Storage Temperature | Expiration Date |
| Proinflammatory Panel 1 (human) Control 1               | A00C0247   | 2-8 $^{\circ}$ C    | 31 MAY 2019     |
| Proinflammatory Panel 1 (human) Control 2               | A00C0248   | 2-8 $^{\circ}$ C    | 31 MAY 2019     |
| Proinflammatory Panel 1 (human) Control 3               | A00C0249   | 2-8 $^{\circ}$ C    | 31 MAY 2019     |

**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> | Dana Short | Quality Control |           | 28 MAY 2014 |