



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Proinflammatory Panel 1 (human) Control Pack

Catalog Number: C4049-1

Lot Number: CP000104

Expiration Date: 31 MAY 2025

Concentration Assignment:

| Analyte | Control 1 | Control 2 | Control 3 | Unit of Measure |
|---------------|-----------|-----------|-----------|-----------------|
| IFN- γ | 612 | 164 | 44.2 | pg/mL |
| IL-1 β | 235 | 55.9 | 11.9 | pg/mL |
| IL-2 | 390 | 39.8 | 3.46 | pg/mL |
| IL-4 | 94.3 | 20.0 | 2.80 | pg/mL |
| IL-6 | 255 | 47.2 | 7.83 | pg/mL |
| IL-8 | 233 | 36.8 | 6.29 | pg/mL |
| IL-10 | 134 | 25.5 | 3.93 | pg/mL |
| IL-12p70 | 177 | 31.9 | 6.39 | pg/mL |
| IL-13 | 164 | 45.2 | 12.1 | pg/mL |
| TNF- α | 96.2 | 16.9 | 2.32 | pg/mL |

Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|-------------------------------------------|------------|---------------------|-----------------|
| Proinflammatory Panel 1 (human) Control 1 | A00C0714 | 2-8°C | 31 MAY 2025 |
| Proinflammatory Panel 1 (human) Control 2 | A00C0715 | 2-8°C | 31 MAY 2025 |
| Proinflammatory Panel 1 (human) Control 3 | A00C0716 | 2-8°C | 30 JUN 2025 |

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|-----------------|--------------------|----------|-----------|-------------|
| Review/Approval | Carissa B. Hubbard | Quality | | 10 JUL 2020 |