



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Cytokine Panel 2 (human) Control Pack

Catalog Number: C4084-1

Lot Number: CP000086

Expiration Date: 31 JAN 2024

Concentration Assignment:

| Analyte | Control 1 | Control 2 | Control 3 | Unit of Measure |
|----------|-----------|-----------|-----------|-----------------|
| IL-17A/F | 5300 | 609 | 78.6 | pg/mL |
| IL-17B | 1480 | 288 | 129 | pg/mL |
| IL-17C | 2120 | 155 | 79.1 | pg/mL |
| IL-17D | 5880 | 512 | 131 | pg/mL |
| IL-1RA | 503 | 88.5 | 51.2 | pg/mL |
| IL-3 | 2870 | 505 | 206 | pg/mL |
| IL-9 | 1290 | 298 | 74.2 | pg/mL |
| TSLP | 618 | 70.9 | 12.7 | pg/mL |

Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|------------------------------------|------------|---------------------|-----------------|
| Cytokine Panel 2 (human) Control 1 | A00C0637 | 2-8°C | 31 JAN 2024 |
| Cytokine Panel 2 (human) Control 2 | A00C0638 | 2-8°C | 31 JAN 2024 |
| Cytokine Panel 2 (human) Control 3 | A00C0639 | 2-8°C | 31 JAN 2024 |

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

| | Name | Function | Signature | Date |
|------------------------|--------------------|----------|-----------|-------------|
| Review/Approval | Carissa B. Hubbard | Quality | | 29 DEC 2020 |