



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** TH17 Panel 1 (mouse) Control Pack

**Catalog Number:** C4246-1

**Lot Number:** CP000062

**Expiration Date:** 30 SEP 2022

**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

| Units = pg/mL           | Assay                |             |             |              |             |             |              |             |              |                    |
|-------------------------|----------------------|-------------|-------------|--------------|-------------|-------------|--------------|-------------|--------------|--------------------|
|                         | ms<br>MIP-3 $\alpha$ | ms<br>IL-22 | ms<br>IL-23 | ms<br>IL-17C | ms<br>IL-31 | ms<br>IL-21 | ms<br>IL-17F | ms<br>IL-16 | ms<br>IL-17A | ms<br>IL-17E/IL-25 |
| Control 1<br>(A00C0522) | 823                  | 445         | 6920        | 1290         | 24 100      | 11 800      | 59 000       | 1760        | 494          | 3860               |
| Control 2<br>(A00C0544) | 200                  | 133         | 1330        | 203          | 4990        | 2100        | 10 500       | 1010        | 85.2         | 1060               |
| Control 3<br>(A00C0543) | 54.6                 | 23.2        | 199         | 35.8         | 1140        | 355         | 2360         | 548         | 14.8         | 283                |

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

| Control Pack Components        |            |                     |                 |
|--------------------------------|------------|---------------------|-----------------|
| Description                    | Lot Number | Storage Temperature | Expiration Date |
| TH17 Panel 1 (mouse) Control 1 | A00C0522   | 2-8°C               | 30 SEP 2022     |
| TH17 Panel 1 (mouse) Control 2 | A00C0544   | 2-8°C               | 30 NOV 2022     |
| TH17 Panel 1 (mouse) Control 3 | A00C0543   | 2-8°C               | 31 OCT 2022     |

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

|                        | Name         | Function          | Signature           | Date        |
|------------------------|--------------|-------------------|---------------------|-------------|
| <b>Review/Approval</b> | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 12 DEC 2017 |