



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Control Pack Description:** TH17 Panel 1 (human) Control Pack

**Catalog Number:** C4085-1

**Lot Number:** CP000090

**Expiration Date:** 30 APR 2024

### Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IL-17A	1980	293	42.5	pg/mL
IL-21	872	161	26.8	pg/mL
IL-31	1080	176	23.9	pg/mL
IL-27	31000	6560	1380	pg/mL
IL-23	5360	602	69.8	pg/mL
IL-22	526	83.3	15.4	pg/mL
MIP-3 $\alpha$	428	175	64.6	pg/mL

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
TH17 Panel 1 (human) Control 1	A00C0657	2-8°C	30 APR 2024
TH17 Panel 1 (human) Control 2	A00C0658	2-8°C	30 APR 2024
TH17 Panel 1 (human) Control 3	A00C0659	2-8°C	30 APR 2024

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		30 MAR 2020