



### CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C4085-1

**Lot Number:** CP000059

**Expiration Date:** 28 FEB 2022

**Storage Temperature:** 2-8°C

**Concentration of Materials Tested:**

Units = pg/mL	Assay						
	hu IL-17A	hu IL-21	hu IL-31	hu IL-27	hu IL-23	hu IL-22	hu MIP-3α
Control 1 (A00C0486)	2690	929	931	18 400	4190	547	531
Control 2 (A00C0477)	429	145	145	3450	487	103	214
Control 3 (A00C0478)	61.4	25.0	21.9	790	58.2	17.7	82.3

**Reconstitution Volume:** 250 µL

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
TH17 Panel 1 (human) Control 1	A00C0486	2-8°C	28 FEB 2022
TH17 Panel 1 (human) Control 2	A00C0477	2-8°C	28 FEB 2022
TH17 Panel 1 (human) Control 3	A00C0478	2-8°C	28 FEB 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>	11 OCT 2018