



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: TH17 Panel 1 (mouse) Control Pack

Catalog Number: C4246-1

Lot Number: CP000060

Expiration Date: 30 SEP 2022

Storage Temperature: 2-8°C

Concentration of Materials Tested:

Units = pg/mL	Assay									
	ms MIP-3α	ms IL-22	ms IL-23	ms IL-17C	ms IL-31	ms IL-21	ms IL-17F	ms IL-16	ms IL-17A	ms IL-17E/IL-25
Control 1 (A00C0534)	797	423	6450	1300	24 500	11 200	51 900	1810	508	4230
Control 2 (A00C0523)	188	134	1270	226	4740	1970	10 800	988	101	1240
Control 3 (A00C0524)	55.3	41.7	272	41.6	1010	400	2370	533	14.6	280

Reconstitution Volume: 250 µL

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
TH17 Panel 1 (mouse) Control 1	A00C0534	2-8°C	30 SEP 2022
TH17 Panel 1 (mouse) Control 2	A00C0523	2-8°C	30 SEP 2022
TH17 Panel 1 (mouse) Control 3	A00C0524	2-8°C	30 SEP 2022

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	04 DEC 2017