



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07622

Product Description: TH17 Panel 1 (human) Control Pack
Catalog Number: C4085-1
Lot Number: CP000043
Expiration Date: 30 SEP 2021

Concentration of Materials Tested:

Units = pg/mL	Assay									
	hu IL-17A	hu IL-17E/IL-25	hu IL-17F	hu IL-21	hu IL-31	hu IL-27	hu IL-23	hu IL-22	hu IL-33	hu MIP-3 α
Reconstituted Control 1 (A00C0436)	1,457	1,250	24,367	590	518	8,470	2,637	279	1,023	245.6
Reconstituted Control 2 (A00C0438)	181	172	4823	82.9	69.6	1960	357	47.7	125	90.4
Reconstituted Control 3 (A00C0437)	28.8	20.2	828	14.5	12.0	477	32.8	9.33	25.7	33.5

Reconstitution Volume: 250 μ L

Control concentrations were assigned across three kit lots.

Proinflammatory Panel 1 (human) Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
TH17 Panel 1 (human) Control 1	A00C0436	2-8°C	30 SEP 2021
TH17 Panel 1 (human) Control 2	A00C0438	2-8°C	30 SEP 2021
TH17 Panel 1 (human) Control 3	A00C0437	2-8°C	30 SEP 2021

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD’s specifications. 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>	23 JAN 2017