



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07891

Product Description: TH17 Panel 1 (human) Control Pack
Catalog Number: C4085-1
Lot Number: CP000044
Expiration Date: 30 NOV 2021

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-3 α
Reconstituted Control 1 (A00C0451)	2 725	886	954	15 017	4 173	537	592
Reconstituted Control 2 (A00C0452)	413	172	148	3 221	464	98	235
Reconstituted Control 3 (A00C0453)	57	26	22	1 055	53	16	83

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	A00C0451	2-8°C	30 NOV 2021
TH17 Panel 1 (human) Control 2	A00C0452	2-8°C	30 NOV 2021
TH17 Panel 1 (human) Control 3	A00C0453	2-8°C	30 NOV 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>	02 MAR 2017