

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Cytokine Panel 2 (human) Control Pack
Catalog Number: C4084-1
Lot Number: CP000180
Expiration Date: 30 JUN 2030

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 2 (human) Control 1	A00C1013	2-8°C	30 JUN 2030
Cytokine Panel 2 (human) Control 2	A00C1014	2-8°C	30 JUN 2030
Cytokine Panel 2 (human) Control 3	A00C1015	2-8°C	30 JUN 2030

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IL-17A/F	4870	600	86.7	pg/mL
IL-17B	2420	459	132	pg/mL
IL-17C	1980	322	110	pg/mL
IL-17D	5930	561	130	pg/mL
IL-1RA	894	157	65.8	pg/mL
IL-3	2920	580	255	pg/mL
IL-9	1310	296	76.8	pg/mL
TSLP	588	80.3	13.1	pg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		04 AUG2025