



### CERTIFICATE OF ANALYSIS

**Control Pack Description:** Cytokine Panel 2 (human) Control Pack  
**Catalog Number:** C4084-1  
**Lot Number:** CP000138  
**Expiration Date:** 31 AUG 2027

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 2 (human) Control 1	A00C0860	2-8°C	31 AUG 2027
Cytokine Panel 2 (human) Control 2	A00C0861	2-8°C	31 AUG 2027
Cytokine Panel 2 (human) Control 3	A00C0862	2-8°C	31 AUG 2027

**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IL-17A/F	5850	718	93.1	pg/mL
IL-17B	3970	373	78.1	pg/mL
IL-17C	2180	337	90.4	pg/mL
IL-17D	7610	811	184	pg/mL
IL-1RA	1010	172	69.4	pg/mL
IL-3	3010	608	257	pg/mL
IL-9	1410	339	84.6	pg/mL
TSLP	668	83.2	14.0	pg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		10 NOV 2023