



### CERTIFICATE OF ANALYSIS

**Control Pack Description:** Cytokine Panel 1 (human) Control Pack  
**Catalog Number:** C4050-1  
**Lot Number:** CP000155  
**Expiration Date:** 30 JUN 2027

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (human) Control 1	A00C0837	2-8°C	30 JUN 2027
Cytokine Panel 1 (human) Control 2	A00C0838	2-8°C	30 JUN 2027
Cytokine Panel 1 (human) Control 3	A00C0920	2-8°C	31 AUG 2028

**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
GM-CSF	456	71.8	7.93	pg/mL
IL-1 $\alpha$	167	58.8	16.8	pg/mL
IL-5	340	117	32.8	pg/mL
IL-7	363	65.1	8.67	pg/mL
IL-12/IL-23p40	951	210	36.1	pg/mL
IL-15	297	48.8	8.57	pg/mL
IL-16	1020	318	98.4	pg/mL
IL-17A	2360	335	54.3	pg/mL
TNF- $\beta$	279	43.6	6.15	pg/mL
VEGF-A	431	144	56.7	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		06 MAR 2024