

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Cytokine Panel 1 (human) Control Pack
Catalog Number: C4050-1
Lot Number: CP000176
Expiration Date: 28 FEB 2030

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (human) Control 1	A00C0977	2-8°C	28 FEB 2030
Cytokine Panel 1 (human) Control 2	A00C0978	2-8°C	28 FEB 2030
Cytokine Panel 1 (human) Control 3	A00C0979	2-8°C	28 FEB 2030

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
GM-CSF	382	47.0	9.22	pg/mL
IL-1 α	146	48.4	13.5	pg/mL
IL-5	335	106	32.1	pg/mL
IL-7	392	61.2	8.68	pg/mL
IL-12/IL-23p40	1360	203	31.7	pg/mL
IL-15	293	42.7	8.11	pg/mL
IL-16	1020	299	85.7	pg/mL
IL-17A	2840	352	50.6	pg/mL
TNF- β	250	42.1	5.41	pg/mL
VEGF-A	390	126	44.4	pg/mL

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

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		25 APR 2025