

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Cytokine Panel 1 (human) Control Pack
Catalog Number: C4050-1
Lot Number: CP000151
Expiration Date: 31 AUG 2028

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (human) Control 1	A00C0918	2-8°C	31 AUG 2028
Cytokine Panel 1 (human) Control 2	A00C0919	2-8°C	31 AUG 2028
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	404	62.0	7.93	pg/mL
IL-1 α	164	54.4	16.8	pg/mL
IL-5	335	109	32.8	pg/mL
IL-7	396	64.5	8.67	pg/mL
IL-12/IL-23p40	1350	215	36.1	pg/mL
IL-15	292	47.5	8.57	pg/mL
IL-16	1080	333	98.4	pg/mL
IL-17A	2800	408	54.3	pg/mL
TNF- β	260	41.2	6.15	pg/mL
VEGF-A	572	190	56.7	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		17 OCT 2023