



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Cytokine Panel 1 (human) Control Pack

Catalog Number: C4050-1

Lot Number: CP000135

Expiration Date: 30 JUN 2027

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
GM-CSF	456	71.8	11.0	pg/mL
IL-1 α	167	58.8	16.3	pg/mL
IL-5	340	117	34.7	pg/mL
IL-7	363	65.1	8.99	pg/mL
IL-12/23p40	951	210	33.4	pg/mL
IL-15	297	48.8	8.63	pg/mL
IL-16	1020	318	109	pg/mL
IL-17A	2360	335	45.4	pg/mL
TNF- β	279	43.6	6.08	pg/mL
VEGF	431	144	47.3	pg/mL

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	A00C0839	2-8°C	30 JUN 2027

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		12 OCT 2022