



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4050-1

Lot Number: CP000091

Expiration Date: 30 APR 2024

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
GM-CSF	394	61.7	8.11	pg/mL
IL-1 α	174	57.6	17.1	pg/mL
IL-5	326	103	32.2	pg/mL
IL-7	342	52.4	8.16	pg/mL
IL-12/23p40	1 320	190	30.4	pg/mL
IL-15	259	40.7	7.82	pg/mL
IL-16	1 270	349	94.7	pg/mL
IL-17A	2 590	374	58.6	pg/mL
TNF- β	241	38.2	5.97	pg/mL
VEGF	472	143	49.6	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (human) Control 1	A00C0646	2-8°C	30 APR 2024
Cytokine Panel 1 (human) Control 2	A00C0647	2-8°C	30 APR 2024
Cytokine Panel 1 (human) Control 3	A00C0648	2-8°C	31 MAY 2024

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		20 JUN 2019