



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4050-1

Lot Number: CP000101

Expiration Date: 30 NOV 2024

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
GM-CSF	435	69.6	11.2	pg/mL
IL-1 α	124	38.5	13.0	pg/mL
IL-5	334	107	33.1	pg/mL
IL-7	346	53.2	9.30	pg/mL
IL-12/23 p40	1490	225	39.8	pg/mL
IL-15	311	50.5	9.58	pg/mL
IL-16	1310	339	103	pg/mL
IL-17A	2220	305	46.4	pg/mL
TNF- β	213	35.4	9.07	pg/mL
VEGF	437	135	45.8	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (human) Control 1	A00C0679	2-8°C	30 NOV 2024
Cytokine Panel 1 (human) Control 2	A00C0680	2-8°C	30 NOV 2024
Cytokine Panel 1 (human) Control 3	A00C0681	2-8°C	30 NOV 2024

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		18 MAR 2020