

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Proinflammatory Panel 1 (human) Control Pack
Catalog Number: C4049-1
Lot Number: CP000184
Expiration Date: 30 SEP 2030

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C1031	2-8°C	30 SEP 2030
Proinflammatory Panel 1 (human) Control 2	A00C1032	2-8°C	30 SEP 2030
Proinflammatory Panel 1 (human) Control 3	A00C1033	2-8°C	30 SEP 2030

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	450	119	31.8	pg/mL
IL-1 β	201	49.3	11.1	pg/mL
IL-2	385	38.9	3.30	pg/mL
IL-4	100	20.6	3.07	pg/mL
IL-6	263	49.9	8.51	pg/mL
IL-8	219	34.3	6.19	pg/mL
IL-10	131	19.2	2.01	pg/mL
IL-12p70	175	31.6	5.03	pg/mL
IL-13	154	47.9	13.3	pg/mL
TNF- α	85.5	16.6	2.52	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		21 NOV 2025