



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4049-1

Lot Number: CP000124

Expiration Date: 31 JUL 2026

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	556	154	39.9	pg/mL
IL-1 β	225	53.4	11.2	pg/mL
IL-2	392	38.0	3.19	pg/mL
IL-4	110	21.4	3.10	pg/mL
IL-6	271	48.9	8.13	pg/mL
IL-8	235	37.0	6.06	pg/mL
IL-10	134	26.3	3.80	pg/mL
IL-12p70	173	31.3	5.64	pg/mL
IL-13	161	46.0	13.2	pg/mL
TNF- α	76.3	14.1	1.77	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0804	2-8°C	31 JUL 2026
Proinflammatory Panel 1 (human) Control 2	A00C0805	2-8°C	31 AUG 2026
Proinflammatory Panel 1 (human) Control 3	A00C0806	2-8°C	31 AUG 2026

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		03 JAN 2022