

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Proinflammatory Panel 1 (human) Control Pack
Catalog Number: C4049-1
Lot Number: CP000177
Expiration Date: 31 MAR 2030

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C1001	2-8°C	31 MAR 2030
Proinflammatory Panel 1 (human) Control 2	A00C1002	2-8°C	31 MAR 2030
Proinflammatory Panel 1 (human) Control 3	A00C1003	2-8°C	31 MAR 2030

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	498	130	35.2	pg/mL
IL-1 β	193	46.4	9.93	pg/mL
IL-2	355	36.2	3.47	pg/mL
IL-4	96.0	19.6	2.59	pg/mL
IL-6	258	47.8	8.27	pg/mL
IL-8	223	34.9	6.64	pg/mL
IL-10	132	24.1	3.28	pg/mL
IL-12p70	170	30.7	5.94	pg/mL
IL-13	183	51.1	14.2	pg/mL
TNF- α	90.1	14.7	2.14	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		06 JUN 2025