

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Proinflammatory Panel 1 (human) Control Pack
Catalog Number: C4049-1
Lot Number: CP000173
Expiration Date: 30 NOV 2029

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0974	2-8°C	30 NOV 2029
Proinflammatory Panel 1 (human) Control 2	A00C0989	2-8°C	31 JAN 2030
Proinflammatory Panel 1 (human) Control 3	A00C0976	2-8°C	30 NOV 2029

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	491	137	36.3	pg/mL
IL-1 β	201	50.9	10.5	pg/mL
IL-2	364	35.5	3.62	pg/mL
IL-4	104	21.7	2.69	pg/mL
IL-6	266	49.5	8.20	pg/mL
IL-8	222	34.2	6.50	pg/mL
IL-10	136	24.7	3.71	pg/mL
IL-12p70	170	32.9	6.08	pg/mL
IL-13	183	52.3	15.1	pg/mL
TNF- α	84.9	15.7	2.12	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		20 MAR 2025