

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Proinflammatory Panel 1 (human) Control Pack
Catalog Number: C4049-1
Lot Number: CP000148
Expiration Date: 30 JUN 2028

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0909	2-8°C	30 JUN 2028
Proinflammatory Panel 1 (human) Control 2	A00C0910	2-8°C	30 JUN 2028
Proinflammatory Panel 1 (human) Control 3	A00C0911	2-8°C	31 JUL 2028

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	533	146	39.9	pg/mL
IL-1 β	245	60.0	12.1	pg/mL
IL-2	433	43.9	4.08	pg/mL
IL-4	97.0	19.8	3.13	pg/mL
IL-6	277	50.0	8.04	pg/mL
IL-8	236	37.6	6.13	pg/mL
IL-10	139	27.0	3.92	pg/mL
IL-12p70	189	34.6	7.09	pg/mL
IL-13	164	45.5	12.6	pg/mL
TNF- α	91.7	15.1	2.38	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		28 AUG 2023