



CERTIFICATE OF ANALYSIS

Control Pack Description: Proinflammatory Panel 1 (human) Control Pack
Catalog Number: C4049-1
Lot Number: CP000154
Expiration Date: 31 JUL 2027

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	A00C0841	2-8°C	31 JUL 2027
Proinflammatory Panel 1 (human) Control 3	A00C0842	2-8°C	31 JUL 2027

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	533	136	40.3	pg/mL
IL-1 β	245	53.8	10.6	pg/mL
IL-2	433	39.5	3.33	pg/mL
IL-4	97.0	20.9	3.02	pg/mL
IL-6	277	48.3	7.51	pg/mL
IL-8	236	39.2	6.62	pg/mL
IL-10	139	26.4	3.90	pg/mL
IL-12p70	189	32.8	6.42	pg/mL
IL-13	164	43.3	11.8	pg/mL
TNF- α	91.7	16.8	2.03	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		07 MAR 2024