



CERTIFICATE OF ANALYSIS

Control Pack Description: Proinflammatory Panel 1 (human) Control Pack
Catalog Number: C4049-1
Lot Number: CP000165
Expiration Date: 30 JUN 2029

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0940	2-8°C	30 JUN 2029
Proinflammatory Panel 1 (human) Control 2	A00C0941	2-8°C	30 JUN 2029
Proinflammatory Panel 1 (human) Control 3	A00C0942	2-8°C	30 JUN 2029

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	569	137	25.8	pg/mL
IL-1 β	241	57.1	12.0	pg/mL
IL-2	368	37.4	3.16	pg/mL
IL-4	88.8	17.8	2.57	pg/mL
IL-6	285	53.4	8.53	pg/mL
IL-8	178	26.0	12.3	pg/mL
IL-10	137	27.3	3.72	pg/mL
IL-12p70	183	33.3	6.13	pg/mL
IL-13	175	49.8	14.1	pg/mL
TNF- α	93.1	16.3	2.34	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		11 APR 2025