



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	587	141	27.4	pg/mL
IL-1 β	237	56.2	12.3	pg/mL
IL-2	391	40.6	3.45	pg/mL
IL-4	97.5	19.7	2.93	pg/mL
IL-6	286	52.9	8.93	pg/mL
IL-8	177	26.0	12.3	pg/mL
IL-10	136	27.1	3.75	pg/mL
IL-12p70	187	33.8	6.24	pg/mL
IL-13	176	50.1	14.6	pg/mL
TNF- α	89.9	15.2	2.23	pg/mL

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Xiomara Talavera	Quality		04 SEP 2024