



CERTIFICATE OF ANALYSIS

Control Pack Description: Proinflammatory Panel 1 (human) Control Pack
Catalog Number: C4049-1
Lot Number: CP000156
Expiration Date: 31 DEC 2028

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0924	2-8°C	31 DEC 2028
Proinflammatory Panel 1 (human) Control 2	A00C0925	2-8°C	31 DEC 2028
Proinflammatory Panel 1 (human) Control 3	A00C0934	2-8°C	31 JAN 2029

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	489	150	35.1	pg/mL
IL-1 β	230	54.9	12.2	pg/mL
IL-2	383	38.3	2.89	pg/mL
IL-4	94.0	18.2	2.27	pg/mL
IL-6	266	51.0	8.11	pg/mL
IL-8	238	36.4	6.68	pg/mL
IL-10	135	27.5	2.81	pg/mL
IL-12p70	172	33.3	5.84	pg/mL
IL-13	183	48.5	13.1	pg/mL
TNF- α	88.1	16.8	1.61	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		29 JUL 2024