



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Proinflammatory Panel 1 (human) Control Pack

Catalog Number: C4049-1

Lot Number: CP000143

Expiration Date: 29 FEB 2028

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	523	139	39.7	pg/mL
IL-1 β	225	55.5	12.2	pg/mL
IL-2	329	38.4	4.40	pg/mL
IL-4	112	20.4	3.02	pg/mL
IL-6	229	47.6	7.82	pg/mL
IL-8	297	35.5	6.09	pg/mL
IL-10	137	24.6	3.85	pg/mL
IL-12p70	177	32.8	5.94	pg/mL
IL-13	156	42.2	12.6	pg/mL
TNF- α	81.5	18.4	2.35	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0870	2-8°C	29 FEB 2028
Proinflammatory Panel 1 (human) Control 2	A00C0871	2-8°C	29 FEB 2028
Proinflammatory Panel 1 (human) Control 3	A00C0872	2-8°C	29 FEB 2028

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 APR 2023