



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4049-1

Lot Number: CP000120

Expiration Date: 31 MAY 2026

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	608	156	41.5	pg/mL
IL-1 β	218	53.0	11.2	pg/mL
IL-2	375	39.0	3.64	pg/mL
IL-4	95.4	19.3	2.71	pg/mL
IL-6	258	47.7	7.49	pg/mL
IL-8	239	38.1	6.47	pg/mL
IL-10	136	26.2	3.77	pg/mL
IL-12p70	169	29.7	5.60	pg/mL
IL-13	161	43.8	12.3	pg/mL
TNF- α	83.0	15.4	1.88	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0782	2-8°C	31 MAY 2026
Proinflammatory Panel 1 (human) Control 2	A00C0783	2-8°C	31 MAY 2026
Proinflammatory Panel 1 (human) Control 3	A00C0784	2-8°C	31 MAY 2026

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		13 JUL 2021