



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4049-1

Lot Number: CP000133

Expiration Date: 31 JAN 2027

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	496	138	30.0	pg/mL
IL-1 β	222	54.0	10.8	pg/mL
IL-2	370	37.2	2.89	pg/mL
IL-4	89.3	19.1	2.48	pg/mL
IL-6	264	46.8	7.30	pg/mL
IL-8	228	36.2	5.39	pg/mL
IL-10	133	26.8	3.47	pg/mL
IL-12p70	131	28.3	7.70	pg/mL
IL-13	148	45.2	11.7	pg/mL
TNF- α	90.3	15.1	1.85	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0818	2-8°C	31 JAN 2027
Proinflammatory Panel 1 (human) Control 2	A00C0819	2-8°C	28 FEB 2027
Proinflammatory Panel 1 (human) Control 3	A00C0820	2-8°C	28 FEB 2027

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 AUG 2022