



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4049-1

Lot Number: CP000084

Expiration Date: 31 JAN 2024

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	530	145	35.3	pg/mL
IL-1 β	209	49.9	10.9	pg/mL
IL-2	409	39.4	3.49	pg/mL
IL-4	104	20.6	3.01	pg/mL
IL-6	267	47.7	7.49	pg/mL
IL-8	218	33.0	6.08	pg/mL
IL-10	132	24.8	3.62	pg/mL
IL-12p70	197	38.8	7.11	pg/mL
IL-13	167	45.7	11.8	pg/mL
TNF- α	87.8	15.7	2.02	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0640	2-8°C	31 JAN 2024
Proinflammatory Panel 1 (human) Control 2	A00C0641	2-8°C	29 FEB 2024
Proinflammatory Panel 1 (human) Control 3	A00C0642	2-8°C	29 FEB 2024

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		14 NOV 2019