



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4049-1

Lot Number: CP000097

Expiration Date: 31 OCT 2024

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	546	167	32.2	pg/mL
IL-1 β	209	49.7	10.0	pg/mL
IL-2	384	37.7	2.96	pg/mL
IL-4	109	20.2	3.11	pg/mL
IL-6	289	49.1	7.89	pg/mL
IL-8	220	33.2	5.42	pg/mL
IL-10	135	24.8	3.62	pg/mL
IL-12p70	158	31.4	5.92	pg/mL
IL-13	121	40.3	13.2	pg/mL
TNF- α	110	16.4	2.37	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0676	2-8°C	31 OCT 2024
Proinflammatory Panel 1 (human) Control 2	A00C0677	2-8°C	31 OCT 2024
Proinflammatory Panel 1 (human) Control 3	A00C0678	2-8°C	31 OCT 2024

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 MAR 2020