



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4048-1

Lot Number: CP000109

Expiration Date: 30 SEP 2025

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	404	30.8	2.25	pg/mL
IL-1 β	629	46.1	4.37	pg/mL
IL-2	825	64.2	5.40	pg/mL
IL-4	434	74.6	14.4	pg/mL
IL-5	376	58.1	9.79	pg/mL
IL-6	1740	192	27.0	pg/mL
KC/GRO	663	107	17.7	pg/mL
IL-10	1010	334	99.4	pg/mL
IL-12p70	14300	3340	828	pg/mL
TNF- α	149	29.6	5.58	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (mouse) Control 1	A00C0736	2-8°C	30 SEP 2025
Proinflammatory Panel 1 (mouse) Control 2	A00C0737	2-8°C	30 SEP 2025
Proinflammatory Panel 1 (mouse) Control 3	A00C0738	2-8°C	31 OCT 2025

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

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