



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4044-1

Lot Number: CP000126

Expiration Date: 31 OCT 2026

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	4610	1450	480	pg/mL
IL-1 β	11200	2900	884	pg/mL
IL-4	834	233	68.8	pg/mL
IL-5	3810	1970	995	pg/mL
IL-6	10000	2910	1010	pg/mL
KC/GRO	811	346	174	pg/mL
IL-10	23300	4950	1430	pg/mL
IL-13	1400	402	123	pg/mL
TNF- α	939	296	90.3	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (rat) Control 1	A00C0808	2-8°C	31 OCT 2026
Proinflammatory Panel 1 (rat) Control 2	A00C0809	2-8°C	31 OCT 2026
Proinflammatory Panel 1 (rat) Control 3	A00C0810	2-8°C	31 OCT 2026

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		14 FEB 2022