



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4049-1

Lot Number: CP000112

Expiration Date: 31 DEC 2025

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	4650	1440	474	pg/mL
IL-1 β	8280	2320	731	pg/mL
IL-4	943	270	81.2	pg/mL
IL-5	3900	1500	749	pg/mL
IL-6	10300	3140	1050	pg/mL
KC/GRO	842	425	210	pg/mL
IL-10	24100	7250	2060	pg/mL
IL-13	1490	411	123	pg/mL
TNF- α	1080	328	104	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (rat) Control 1	A00C0762	2-8°C	31 DEC 2025
Proinflammatory Panel 1 (rat) Control 2	A00C0763	2-8°C	31 DEC 2025
Proinflammatory Panel 1 (rat) Control 3	A00C0764	2-8°C	31 DEC 2025

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		23 FEB 2021