



CERTIFICATE OF ANALYSIS

Control Pack Description: Proinflammatory Panel 1 (rat) Control Pack
Catalog Number: C4044-1
Lot Number: CP000181
Expiration Date: 30 APR 2030

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (rat) Control 1	A00C0983	2-8°C	30 APR 2030
Proinflammatory Panel 1 (rat) Control 2	A00C0984	2-8°C	30 APR 2030
Proinflammatory Panel 1 (rat) Control 3	A00C0985	2-8°C	30 APR 2030

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	4580	1460	449	pg/mL
IL-1 β	25300	6230	1870	pg/mL
IL-4	1030	296	92.7	pg/mL
IL-5	3580	1850	952	pg/mL
IL-6	11700	3660	1190	pg/mL
KC/GRO	931	469	250	pg/mL
IL-10	16000	4690	1480	pg/mL
IL-13	1330	397	144	pg/mL
TNF- α	900	288	92.7	pg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 AUG 2025