



CERTIFICATE OF ANALYSIS

Control Pack Description: Proinflammatory Panel 1 (rat) Control Pack
Catalog Number: C4044-1
Lot Number: CP000161
Expiration Date: 31 MAY 2029

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (rat) Control 1	A00C0950	2-8°C	31 MAY 2029
Proinflammatory Panel 1 (rat) Control 2	A00C0951	2-8°C	31 MAY 2029
Proinflammatory Panel 1 (rat) Control 3	A00C0952	2-8°C	31 MAY 2029

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	4380	1400	401	pg/mL
IL-1 β	9360	2810	894	pg/mL
IL-4	1040	312	98.3	pg/mL
IL-5	4420	2370	1200	pg/mL
IL-6	12700	4060	1260	pg/mL
KC/GRO	955	475	238	pg/mL
IL-10	18400	5610	1490	pg/mL
IL-13	1480	438	137	pg/mL
TNF- α	943	300	97.3	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		09 JUL 2024