



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4044-1

Lot Number: CP000100

Expiration Date: 30 NOV 2024

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	4580	1450	470	pg/mL
IL-1 β	13200	4150	1420	pg/mL
IL-4	926	262	80.9	pg/mL
IL-5	3550	1740	889	pg/mL
IL-6	9830	3150	1100	pg/mL
KC/GRO	882	403	211	pg/mL
IL-10	25000	7380	1910	pg/mL
IL-13	1430	413	132	pg/mL
TNF- α	948	293	113	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (rat) Control 1	A00C0685	2-8°C	30 NOV 2024
Proinflammatory Panel 1 (rat) Control 2	A00C0686	2-8°C	30 NOV 2024
Proinflammatory Panel 1 (rat) Control 3	A00C0687	2-8°C	30 NOV 2024

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		31 JAN 2020