



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4294-1

Lot Number: CP000131

Expiration Date: 31 MAR 2027

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	119	21.5	1.46	pg/mL
IL-1 β	2370	255	79.8	pg/mL
IL-2	240	40.8	6.50	pg/mL
IL-4	656	110	11.6	pg/mL
IL-5	276	46.4	11.7	pg/mL
IL-6	44400	8110	1900	pg/mL
KC/GRO	645	120	22.0	pg/mL
IL-10	2090	382	73.3	pg/mL
IL-13	1520	442	130	pg/mL
TNF- α	859	268	90.4	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (rat) Control 1	A00C0815	2-8°C	31 MAR 2027
Proinflammatory Panel 1 (rat) Control 2	A00C0816	2-8°C	31 MAR 2027
Proinflammatory Panel 1 (rat) Control 3	A00C0817	2-8°C	31 MAR 2027

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		09 MAY 2022