



CERTIFICATE OF ANALYSIS

Control Pack Description: Proinflammatory Panel 1 (rat) Control Pack
Catalog Number: C4294-1
Lot Number: CP000157
Expiration Date: 31 MAR 2029

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (rat) Control 1	A00C0927	2-8°C	31 MAR 2029
Proinflammatory Panel 1 (rat) Control 2	A00C0928	2-8°C	31 MAR 2029
Proinflammatory Panel 1 (rat) Control 3	A00C0929	2-8°C	31 MAR 2029

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	390	75.2	14.9	pg/mL
IL-1 β	865	99.7	14.8	pg/mL
IL-2	639	96.3	14.5	pg/mL
IL-4	519	156	44.2	pg/mL
IL-5	460	113	24.4	pg/mL
IL-6	73700	11300	2020	pg/mL
KC/GRO	862	136	18.9	pg/mL
IL-10	6140	1980	591	pg/mL
IL-13	1480	464	159	pg/mL
TNF- α	1140	136	17.1	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		29 APR 2024