



CERTIFICATE OF ANALYSIS

Control Pack Description: Proinflammatory Panel 1 (rat) Control Pack
Catalog Number: C4044-1
Lot Number: CP000142
Expiration Date: 31 JAN 2028

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	4420	1500	487	pg/mL
IL-1 β	8730	3680	1400	pg/mL
IL-4	866	244	78.1	pg/mL
IL-5	3610	1870	990	pg/mL
IL-6	9680	3200	887	pg/mL
KC/GRO	762	366	183	pg/mL
IL-10	16800	5380	1580	pg/mL
IL-13	1300	362	114	pg/mL
TNF- α	966	314	95.5	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (rat) Control 1	A00C0867	2-8°C	31 JAN 2028
Proinflammatory Panel 1 (rat) Control 2	A00C0868	2-8°C	31 JAN 2028
Proinflammatory Panel 1 (rat) Control 3	A00C0869	2-8°C	31 JAN 2028

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 MAR 2023