

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Proinflammatory Panel 1 (mouse) Control Pack
Catalog Number: C4048-1
Lot Number: CP000158
Expiration Date: 28 FEB 2029

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (mouse) Control 1	A00C0931	2-8°C	28 FEB 2029
Proinflammatory Panel 1 (mouse) Control 2	A00C0932	2-8°C	28 FEB 2029
Proinflammatory Panel 1 (mouse) Control 3	A00C0933	2-8°C	28 FEB 2029

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- γ	504	34.4	2.36	pg/mL
IL-1 β	782	55.0	4.81	pg/mL
IL-2	951	67.7	5.67	pg/mL
IL-4	586	91.0	16.5	pg/mL
IL-5	398	60.6	9.48	pg/mL
IL-6	1680	251	34.6	pg/mL
KC/GRO	842	110	17.8	pg/mL
IL-10	1060	323	121	pg/mL
IL-12p70	19800	3840	581	pg/mL
TNF- α	113	28.3	5.43	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		13 MAY 2024