



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Control Pack Description:** Proinflammatory Panel 1 (mouse) Control Pack

**Catalog Number:** C4048-1

**Lot Number:** CP000102

**Expiration Date:** 30 APR 2023

### Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IFN- $\gamma$	469	32.0	2.12	pg/mL
IL-1 $\beta$	722	52.0	4.32	pg/mL
IL-2	908	54.5	5.57	pg/mL
IL-4	479	62.1	14.5	pg/mL
IL-5	356	54.2	8.30	pg/mL
IL-6	990	152	28.9	pg/mL
KC/GRO	709	119	17.3	pg/mL
IL-10	1100	331	124	pg/mL
IL-12p70	14400	3710	822	pg/mL
TNF- $\alpha$	158	28.1	4.46	pg/mL

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (mouse) Control 1	A00C0717	2-8°C	30 APR 2023
Proinflammatory Panel 1 (mouse) Control 2	A00C0718	2-8°C	30 APR 2023
Proinflammatory Panel 1 (mouse) Control 3	A00C0719	2-8°C	30 APR 2023

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		27 MAY 2020