



## RFE MEMO


Page 1 of 1

**Product Description(s):** Cytokine Panel 1 (mouse) Control 1

**Catalog Number(s) / Size(s):** 3-3081-420181-B

**Lot Number:** A00C0491

Based on additional testing, the original date of expiration for this control has been extended by 12 months. The expiration date of A00C0491 has been extended to 31 MAR 2023.

Function	Name	Signature	Date
Quality	Karen Hamilla		01 APR 2022

This memo supersedes all other statements regarding date of expiration for this product.



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C4245-1

**Lot Number:** CP000128

**Expiration Date:** 31 MAR 2023


**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IL-9	3569	822	216	pg/mL
MCP-1	458	145	57.9	pg/mL
IL-33	2382	239	32.9	pg/mL
IL-27p28/IL-30	8677	934	140	pg/mL
IL-15	28253	5150	786	pg/mL
IL-17A/F	2185	206	16.3	pg/mL
MIP-1 $\alpha$	573	43.2	3.95	pg/mL
IP-10	887	252	81.6	pg/mL
MIP-2	567	70.8	10.3	pg/mL

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (mouse) Control 1	A00C0491	2-8°C	31 MAR 2023
Cytokine Panel 1 (mouse) Control 2	A00C0797	2-8°C	30 JUN 2026
Cytokine Panel 1 (mouse) Control 3	A00C0798	2-8°C	30 JUN 2026

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		05 APR 2022