



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4245-1

Lot Number: CP000129

Expiration Date: 30 JUN 2026

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IL-9	3650	822	216	pg/mL
MCP-1	514	145	57.9	pg/mL
IL-33	2190	239	32.9	pg/mL
IL-27p28/IL-30	7990	934	140	pg/mL
IL-15	29700	5150	786	pg/mL
IL-17A/F	2240	206	16.3	pg/mL
MIP-1 α	625	43.2	3.95	pg/mL
IP-10	1010	252	81.6	pg/mL
MIP-2	629	70.8	10.3	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (mouse) Control 1	A00C0821	2-8°C	31 JAN 2027
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
Review/Approval	Karen Hamilla	Quality		13 APR 2022