

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Cytokine Panel 1 (mouse) Control Pack
Catalog Number: C4245-1
Lot Number: CP000186
Expiration Date: 31 MAY 2028

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (mouse) Control 1	A00C1039	2-8°C	30 NOV 2030
Cytokine Panel 1 (mouse) Control 2	A00C1040	2-8°C	30 NOV 2030
Cytokine Panel 1 (mouse) Control 3	A00C0907	2-8°C	31 MAY 2028

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IL-9	3480	863	231	pg/mL
MCP-1	459	142	65.2	pg/mL
IL-33	1950	273	36.9	pg/mL
IL-27p28/IL-30	8970	1630	182	pg/mL
IL-15	28700	4650	740	pg/mL
IL-17A/F	2180	246	19.3	pg/mL
MIP-1 α	562	44.6	4.66	pg/mL
IP-10	980	274	84.2	pg/mL
MIP-2	557	72.1	9.75	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		02 FEB 2026