



CERTIFICATE OF ANALYSIS

Control Pack Description: Cytokine Panel 1 (mouse) Control Pack
Catalog Number: C4245-1
Lot Number: CP000164
Expiration Date: 31 MAR 2028

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (mouse) Control 1	A00C0905	2-8°C	31 MAY 2028
Cytokine Panel 1 (mouse) Control 2	A00C0906	2-8°C	31 MAY 2028
Cytokine Panel 1 (mouse) Control 3	A00C0891	2-8°C	31 MAR 2028

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IL-9	3650	852	211	pg/mL
MCP-1	555	162	61.0	pg/mL
IL-33	2170	248	36.2	pg/mL
IL-27p28/IL-30	7370	1140	124	pg/mL
IL-15	28700	4540	757	pg/mL
IL-17A/F	2330	210	19.1	pg/mL
MIP-1 α	652	53.3	4.09	pg/mL
IP-10	1030	274	84.1	pg/mL
MIP-2	583	67.4	9.49	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		30 AUG 2024