

**CERTIFICATE OF ANALYSIS****Control Pack Description:** Cytokine Panel 1 (mouse) Control Pack**Catalog Number:** C4245-1**Lot Number:** CP000145**Expiration Date:** 31 JAN 2027**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IL-9	3650	846	211	pg/mL
MCP-1	514	137	61.0	pg/mL
IL-33	2190	233	36.2	pg/mL
IL-27p28/IL-30	7990	1220	124	pg/mL
IL-15	29700	4550	757	pg/mL
IL-17A/F	2240	210	19.1	pg/mL
MIP-1 α	625	44.5	4.09	pg/mL
IP-10	1010	269	84.1	pg/mL
MIP-2	629	65.5	9.49	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	A00C0889	2-8°C	31 MAR 2028
Cytokine Panel 1 (mouse) Control 3	A00C0891	2-8°C	31 MAR 2028

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		04 MAY 2023