



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C4245-1

Lot Number: CP000073

Expiration Date: 31 MAR 2023

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
IL-9	3480	809	204	pg/mL
MCP-1	439	133	55.3	pg/mL
IL-33	2350	274	35.8	pg/mL
IL-27p28/IL-30	6980	1030	141	pg/mL
IL-15	27400	4210	724	pg/mL
IL-17A/F	2130	205	17.7	pg/mL
MIP-1 α	771	64.2	5.27	pg/mL
IP-10	857	228	69.8	pg/mL
MIP-2	549	65.0	8.81	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (mouse) Control 1	A00C0538	2-8°C	31 MAR 2023
Cytokine Panel 1 (mouse) Control 2	A00C0539	2-8°C	31 MAR 2023
Cytokine Panel 1 (mouse) Control 3	A00C0540	2-8°C	31 MAR 2023

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		26 FEB 2020