



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 1 (mouse) Control Pack

Catalog Number: C4245-1

Lot Number: CP000051

Expiration Date: 31 MAR 2022

Concentration of Materials Tested:

Units = pg/mL	Assay								
	ms IL-9	ms MCP-1	ms IL-33	ms IL-27p28/IL-30	ms IL-15	ms IL-17A/F	ms MIP-1 α	ms IP-10	ms MIP-2
Reconstituted Control 1 (A00C0491)	3569	458	2382	8677	28 253	2185	573	887	567
Reconstituted Control 2 (A00C0492)	872	150	284	1227	4168	203	49.3	229	66.6
Reconstituted Control 3 (A00C0493)	202	53.6	36.2	187	663	17.9	4.24	59.9	10.1

Reconstitution Volume: 250 μ L

Control Pack 1 Components			
Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (mouse) Control 1	A00C0491	2-8°C	31 MAR 2022
Cytokine Panel 1 (mouse) Control 2	A00C0492	2-8°C	31 MAR 2022
Cytokine Panel 1 (mouse) Control 3	A00C0493	2-8°C	31 MAR 2022

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	28 JUL 2017