



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 1 (mouse) Calibrator Blend

Catalog Number: C0245-2

Lot Number: A0080323

Expiration Date: 31 MAR 2027

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-9	3880	pg/mL
MCP-1	545	pg/mL
IL-33	3020	pg/mL
IL-27p28/IL-30	9070	pg/mL
IL-15	39200	pg/mL
IL-17A/F	2560	pg/mL
MIP-1 α	528	pg/mL
IP-10	1030	pg/mL
MIP-2	704	pg/mL

Reconstitution Volume: 1000 μ L

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		09 MAY 2022