



CERTIFICATE OF ANALYSIS

Product Description: Cytokine Panel 1 (mouse) Kit Calibrator
Catalog Number: C0245-2
Lot Number: A0080353
Expiration Date: 29 FEB 2028
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-9	3940	pg/mL
MCP-1	479	pg/mL
IL-33	2860	pg/mL
IL-27p28/IL-30	7500	pg/mL
IL-15	41000	pg/mL
IL-17A/F	2530	pg/mL
MIP-1α	537	pg/mL
IP-10	1010	pg/mL
MIP-2	616	pg/mL

Reconstitution Volume: 1000 µL

This product was manufactured and tested according to MSD documents. This lot meets MSD’s specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		15 MAR 2023