



CERTIFICATE OF ANALYSIS

Product Description: Cytokine Panel 1 (mouse) Kit Calibrator
Catalog Number: C0245-2
Lot Number: A0080371
Expiration Date: 30 JUN 2028
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-9	4160	pg/mL
MCP-1	574	pg/mL
IL-33	3220	pg/mL
IL-27p28/IL-30	9910	pg/mL
IL-15	41900	pg/mL
IL-17A/F	2890	pg/mL
MIP-1 α	630	pg/mL
IP-10	1030	pg/mL
MIP-2	692	pg/mL

Reconstitution Volume: 1000 μ L

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 AUG 2023