



## CERTIFICATE OF ANALYSIS

**Product Description:** Cytokine Panel 1 (mouse) Kit Calibrator  
**Catalog Number:** C0245-2  
**Lot Number:** A0080384  
**Expiration Date:** 28 FEB 2029  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-9	3950	pg/mL
MCP-1	528	pg/mL
IL-33	3010	pg/mL
IL-27p28/IL-30	7150	pg/mL
IL-15	38000	pg/mL
IL-17A/F	2680	pg/mL
MIP-1 $\alpha$	590	pg/mL
IP-10	945	pg/mL
MIP-2	612	pg/mL

**Reconstitution Volume:** 1000  $\mu$ L

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		02 APR 2024