



### CERTIFICATE OF ANALYSIS

**Product Description:** Cytokine Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0050-2  
**Lot Number:** A0080463  
**Expiration Date:** 31 MAR 2031  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GM-CSF	1010	pg/mL
IL-1 $\alpha$	463	pg/mL
IL-5	903	pg/mL
IL-7	853	pg/mL
IL-12/IL-23p40	3590	pg/mL
IL-15	873	pg/mL
IL-16	3550	pg/mL
IL-17A	5020	pg/mL
TNF- $\beta$	703	pg/mL
VEGF-A	1270	pg/mL

**Reconstitution Volume:** 1000  $\mu$ 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
<b>Review/Approval</b>	Karen Hamilla	Quality		12 MAY 2026