

**CERTIFICATE OF ANALYSIS**


**Product Description:** Cytokine Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0050-2  
**Lot Number:** A0080454  
**Expiration Date:** 31 DEC 2030  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GM-CSF	1020	pg/mL
IL-1 $\alpha$	435	pg/mL
IL-5	867	pg/mL
IL-7	944	pg/mL
IL-12/IL-23p40	3530	pg/mL
IL-15	846	pg/mL
IL-16	2550	pg/mL
IL-17A	6310	pg/mL
TNF- $\beta$	752	pg/mL
VEGF-A	1150	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		22 JAN 2026