

**CERTIFICATE OF ANALYSIS**


Product Description: Cytokine Panel 1 (human) Calibrator Blend
Catalog Number: C0050-2
Lot Number: A0080435
Expiration Date: 31 MAY 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GM-CSF	966	pg/mL
IL-1 α	435	pg/mL
IL-5	881	pg/mL
IL-7	878	pg/mL
IL-12/IL-23p40	3570	pg/mL
IL-15	889	pg/mL
IL-16	2560	pg/mL
IL-17A	5800	pg/mL
TNF- β	677	pg/mL
VEGF-A	1120	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		04 AUG 2025