



CERTIFICATE OF ANALYSIS

Product Description: Cytokine Panel 1 (human) Calibrator Blend
Catalog Number: C0050-2
Lot Number: A0080374
Expiration Date: 31 JUL 2028
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GM-CSF	1080	pg/mL
IL-1 α	407	pg/mL
IL-5	881	pg/mL
IL-7	881	pg/mL
IL-12/IL-23p40	3530	pg/mL
IL-15	821	pg/mL
IL-16	2990	pg/mL
IL-17A	7360	pg/mL
TNF- β	687	pg/mL
VEGF-A	1660	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		12 SEP 2023