



CERTIFICATE OF ANALYSIS

**Product Description:** Cytokine Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0050-2  
**Lot Number:** A0080383  
**Expiration Date:** 31 DEC 2028  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GM-CSF	1110	pg/mL
IL-1 $\alpha$	412	pg/mL
IL-5	826	pg/mL
IL-7	887	pg/mL
IL-12/IL-23p40	3460	pg/mL
IL-15	850	pg/mL
IL-16	2360	pg/mL
IL-17A	6490	pg/mL
TNF- $\beta$	721	pg/mL
VEGF-A	1820	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>	Michelle Likiam	Quality		25 JAN 2024