

**CERTIFICATE OF ANALYSIS**


Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080427
Expiration Date: 28 FEB 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1420	pg/mL
IL-1 β	570	pg/mL
IL-2	1290	pg/mL
IL-4	240	pg/mL
IL-6	730	pg/mL
IL-8	534	pg/mL
IL-10	380	pg/mL
IL-12p70	495	pg/mL
IL-13	560	pg/mL
TNF- α	362	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		07 APR 2025