



CERTIFICATE OF ANALYSIS


Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080423
Expiration Date: 28 FEB 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1340	pg/mL
IL-10	365	pg/mL
IL-12p70	459	pg/mL
IL-13	570	pg/mL
IL-1 β	549	pg/mL
IL-2	1260	pg/mL
IL-4	241	pg/mL
IL-6	716	pg/mL
IL-8	538	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		13 MAR 2025