



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080461
Expiration Date: 31 MAR 2031
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1600	pg/mL
IL-1 β	600	pg/mL
IL-2	1720	pg/mL
IL-4	255	pg/mL
IL-6	802	pg/mL
IL-8	581	pg/mL
IL-10	374	pg/mL
IL-12p70	495	pg/mL
IL-13	527	pg/mL
TNF- α	346	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		27 APR 2026