



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080466
Expiration Date: 31 JUL 2031
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1630	pg/mL
IL-1 β	613	pg/mL
IL-2	1450	pg/mL
IL-4	240	pg/mL
IL-6	819	pg/mL
IL-8	599	pg/mL
IL-10	376	pg/mL
IL-12p70	482	pg/mL
IL-13	534	pg/mL
TNF- α	342	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		01 SEP 2026