



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080412
Expiration Date: 31 DEC 2029
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1440	pg/mL
IL-1 β	585	pg/mL
IL-2	1390	pg/mL
IL-4	227	pg/mL
IL-6	734	pg/mL
IL-8	575	pg/mL
IL-10	393	pg/mL
IL-12p70	489	pg/mL
IL-13	572	pg/mL
TNF- α	371	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		21 JAN 2025