



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080464
Expiration Date: 30 APR 2031
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1520	pg/mL
IL-1 β	595	pg/mL
IL-2	1720	pg/mL
IL-4	254	pg/mL
IL-6	796	pg/mL
IL-8	583	pg/mL
IL-10	375	pg/mL
IL-12p70	479	pg/mL
IL-13	531	pg/mL
TNF- α	344	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		22 JUN 2026