

**CERTIFICATE OF ANALYSIS**


Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080434
Expiration Date: 31 MAY 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1170	pg/mL
IL-1 β	622	pg/mL
IL-2	1560	pg/mL
IL-4	276	pg/mL
IL-6	780	pg/mL
IL-8	555	pg/mL
IL-10	381	pg/mL
IL-12p70	494	pg/mL
IL-13	596	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		11 AUG 2025