

**CERTIFICATE OF ANALYSIS**


Product Description: Proinflammatory Panel 1 (rat) Calibrator Blend
Catalog Number: C0044-2
Lot Number: A0080413
Expiration Date: 30 APR 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	6300	pg/mL
IL-1 β	11900	pg/mL
IL-4	1050	pg/mL
IL-5	9540	pg/mL
IL-6	13400	pg/mL
KC/GRO	2950	pg/mL
IL-10	23900	pg/mL
IL-13	1670	pg/mL
TNF- α	1330	pg/mL

Reconstitution Volume: 1000 μ L

This product was manufactured and tested according to MSD documents. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		10 JUN 2025