

**CERTIFICATE OF ANALYSIS**


Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend
Catalog Number: C0048-2
Lot Number: A0080429
Expiration Date: 31 MAY 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	938	pg/mL
IL-1 β	1650	pg/mL
IL-2	2410	pg/mL
IL-4	1810	pg/mL
IL-5	907	pg/mL
IL-6	4680	pg/mL
KC/GRO	2210	pg/mL
IL-10	2870	pg/mL
IL-12p70	29800	pg/mL
TNF- α	630	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		03 JUL 2025