



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend
Catalog Number: C0048-2
Lot Number: A0080351
Expiration Date: 31 DEC 2027
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1030	pg/mL
IL-1 β	1740	pg/mL
IL-2	2580	pg/mL
IL-4	1760	pg/mL
IL-5	1000	pg/mL
IL-6	4930	pg/mL
KC/GRO	2320	pg/mL
IL-10	3860	pg/mL
IL-12p70	27900	pg/mL
TNF- α	622	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		10 MAR 2023