



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (rat) Calibrator Blend
Catalog Number: C0044-2
Lot Number: A0080352
Expiration Date: 31 DEC 2027
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	5287	pg/mL
IL-1 β	22200	pg/mL
IL-4	1133	pg/mL
IL-5	10153	pg/mL
IL-6	13900	pg/mL
KC/GRO	2683	pg/mL
IL-10	25900	pg/mL
IL-13	1693	pg/mL
TNF- α	1283	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		08 MAR 2023