



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (rat) Calibrator Blend
Catalog Number: C0044-2
Lot Number: A0080386
Expiration Date: 28 FEB 2029
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	5230	pg/mL
IL-1 β	12700	pg/mL
IL-4	1110	pg/mL
IL-5	8440	pg/mL
IL-6	13900	pg/mL
KC/GRO	2700	pg/mL
IL-10	30000	pg/mL
IL-13	1560	pg/mL
TNF- α	1200	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 APR 2024