



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (rat) Calibrator Blend
Catalog Number: C0044-2
Lot Number: A0080379
Expiration Date: 30 NOV 2028
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	5430	pg/mL
IL-1 β	13400	pg/mL
IL-4	1140	pg/mL
IL-5	9980	pg/mL
IL-6	15500	pg/mL
KC/GRO	2920	pg/mL
IL-10	29100	pg/mL
IL-13	1630	pg/mL
TNF- α	1220	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 JAN 2024