



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (rat) Calibrator Blend

Catalog Number: C0044-2

Lot Number: A0080256

Expiration Date: 31 MAY 2025

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	5800	pg/mL
IL-1 β	13500	pg/mL
IL-4	1100	pg/mL
IL-5	10800	pg/mL
IL-6	12600	pg/mL
KC/GRO	2650	pg/mL
IL-10	23300	pg/mL
IL-13	1750	pg/mL
TNF- α	1320	pg/mL

Reconstitution Volume: 1000 μ L

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		10 JUN 2020