



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (rat) Calibrator Blend

Catalog Number: C0044-2

Lot Number: A0080308

Expiration Date: 30 SEP 2026

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	6260	pg/mL
IL-1 β	16100	pg/mL
IL-4	1110	pg/mL
IL-5	9230	pg/mL
IL-6	13500	pg/mL
KC/GRO	2570	pg/mL
IL-10	31800	pg/mL
IL-13	1680	pg/mL
TNF- α	1200	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		14 DEC 2021