



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (rat) Calibrator Blend

Catalog Number: C0044-2

Lot Number: A0080337

Expiration Date: 30 SEP 2027

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	5240	pg/mL
IL-1 β	20500	pg/mL
IL-4	988	pg/mL
IL-5	9030	pg/mL
IL-6	13500	pg/mL
KC/GRO	2450	pg/mL
IL-10	27700	pg/mL
IL-13	1500	pg/mL
TNF- α	1150	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		31 OCT 2022