



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend

Catalog Number: C0048-2

Lot Number: A0080329

Expiration Date: 31 MAY 2027

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	890	pg/mL
IL-1 β	1590	pg/mL
IL-2	2260	pg/mL
IL-4	1660	pg/mL
IL-5	947	pg/mL
IL-6	4680	pg/mL
KC/GRO	1780	pg/mL
IL-10	3030	pg/mL
IL-12p70	27500	pg/mL
TNF- α	594	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		03 JUL 2022