



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend

Catalog Number: C0048-2

Lot Number: A0080322

Expiration Date: 28 FEB 2027

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	909	pg/mL
IL-1 β	1765	pg/mL
IL-2	2595	pg/mL
IL-4	1680	pg/mL
IL-5	1027	pg/mL
IL-6	5440	pg/mL
KC/GRO	1960	pg/mL
IL-10	3930	pg/mL
IL-12p70	31400	pg/mL
TNF- α	619	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		30 MAR 2022