



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend

Catalog Number: C0048-2

Lot Number: A0080252

Expiration Date: 30 APR 2025

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	785	pg/mL
IL-1 β	1530	pg/mL
IL-2	2670	pg/mL
IL-4	1600	pg/mL
IL-5	991	pg/mL
KC/GRO	2030	pg/mL
IL-10	2530	pg/mL
IL-12p70	31500	pg/mL
TNF- α	516	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		08 SEP 2020