



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080314

Expiration Date: 31 DEC 2026

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1410	pg/ml
IL-1 β	588	pg/ml
IL-2	1380	pg/ml
IL-4	220	pg/ml
IL-6	715	pg/ml
IL-8	597	pg/ml
IL-10	383	pg/ml
IL-12p70	483	pg/ml
IL-13	456	pg/ml
TNF- α	351	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 FEB 2022