



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080319

Expiration Date: 28 FEB 2027

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1390	pg/mL
IL-1 β	627	pg/mL
IL-2	1550	pg/mL
IL-4	233	pg/mL
IL-6	720	pg/mL
IL-8	615	pg/mL
IL-10	374	pg/mL
IL-12p70	440	pg/mL
IL-13	493	pg/mL
TNF- α	395	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		06 APR 2022