



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080281

Expiration Date: 28 FEB 2026


Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1580	pg/mL
IL-1 β	609	pg/mL
IL-2	1500	pg/mL
IL-4	238	pg/mL
IL-6	737	pg/mL
IL-8	626	pg/mL
IL-10	393	pg/mL
IL-12p70	438	pg/mL
IL-13	482	pg/mL
TNF- α	349	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	Carissa B. Hubbard	Quality		06 APR 2021