



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080305

Expiration Date: 30 SEP 2026

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1340	pg/mL
IL-1 β	613	pg/mL
IL-2	1440	pg/mL
IL-4	266	pg/mL
IL-6	739	pg/mL
IL-8	658	pg/mL
IL-10	384	pg/mL
IL-12p70	480	pg/mL
IL-13	503	pg/mL
TNF- α	337	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		18 NOV 2021