



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080304

Expiration Date: 31 AUG 2026

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1510	pg/mL
IL-1 β	618	pg/mL
IL-2	1410	pg/mL
IL-4	251	pg/mL
IL-6	708	pg/mL
IL-8	619	pg/mL
IL-10	391	pg/mL
IL-12p70	474	pg/mL
IL-13	486	pg/mL
TNF- α	334	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 OCT 2021