



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080267

Expiration Date: 31 AUG 2025

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1360	pg/mL
IL-1 β	601	pg/mL
IL-2	1440	pg/mL
IL-4	227	pg/mL
IL-6	706	pg/mL
IL-8	566	pg/mL
IL-10	377	pg/mL
IL-12p70	474	pg/mL
IL-13	492	pg/mL
TNF- α	397	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		24 AUG 2020