



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080296

Expiration Date: 31 MAY 2026

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1450	pg/mL
IL-1 β	584	pg/mL
IL-2	1390	pg/mL
IL-4	208	pg/mL
IL-6	683	pg/mL
IL-8	580	pg/mL
IL-10	381	pg/mL
IL-12p70	447	pg/mL
IL-13	481	pg/mL
TNF- α	324	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		23 JUL 2021