



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Proinflammatory Panel 1 (human) Calibrator Blend

**Catalog Number:** C0049-2

**Lot Number:** A0080289

**Expiration Date:** 31 MAR 2026

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	1530	pg/mL
IL-1 $\beta$	622	pg/mL
IL-2	1500	pg/mL
IL-4	225	pg/mL
IL-6	756	pg/mL
IL-8	617	pg/mL
IL-10	391	pg/mL
IL-12p70	464	pg/mL
IL-13	493	pg/mL
TNF- $\alpha$	337	pg/mL

**Reconstitution Volume:** 1000  $\mu$ L

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		02 JUN 2021