



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0049-2

**Lot Number:** A0080299

**Expiration Date:** 30 JUN 2026

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	1500	pg/mL
IL-1 $\beta$	601	pg/mL
IL-2	1440	pg/mL
IL-4	242	pg/mL
IL-6	703	pg/mL
IL-8	601	pg/mL
IL-10	386	pg/mL
IL-12p70	472	pg/mL
IL-13	475	pg/mL
TNF- $\alpha$	330	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
<b>Review/Approval</b>	Karen Hamilla	Quality		15 SEP 2021