



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0049-2

**Lot Number:** A0080333

**Expiration Date:** 30 JUN 2027

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	1170	pg/mL
IL-1 $\beta$	559	pg/mL
IL-2	1470	pg/mL
IL-4	232	pg/mL
IL-6	684	pg/mL
IL-8	557	pg/mL
IL-10	343	pg/mL
IL-12p70	427	pg/mL
IL-13	470	pg/mL
TNF- $\alpha$	349	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		12 AUG 2022