



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0048-2

**Lot Number:** A0080264

**Expiration Date:** 31 JUL 2025

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	764	pg/mL
IL-1 $\beta$	1530	pg/mL
IL-2	2530	pg/mL
IL-4	1790	pg/mL
IL-5	994	pg/mL
IL-6	5240	pg/mL
KC/GRO	2020	pg/mL
IL-10	3540	pg/mL
IL-12p70	32500	pg/mL
TNF- $\alpha$	639	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		08 SEP 2020