



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0048-2

**Lot Number:** A0080250

**Expiration Date:** 28 FEB 2025

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	764	pg/mL
IL-1 $\beta$	1510	pg/mL
IL-2	2430	pg/mL
IL-4	1660	pg/mL
IL-5	945	pg/mL
IL-6	5120	pg/mL
KC/GRO	2150	pg/mL
IL-10	2630	pg/mL
IL-12p70	29800	pg/mL
TNF- $\alpha$	572	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		05 MAR 2020