



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0048-2

**Lot Number:** A0080282

**Expiration Date:** 31 MAR 2026

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	742	pg/mL
IL-1 $\beta$	1810	pg/mL
IL-2	2500	pg/mL
IL-4	1660	pg/mL
IL-5	985	pg/mL
IL-6	5070	pg/mL
KC/GRO	1940	pg/mL
IL-10	3240	pg/mL
IL-12p70	31600	pg/mL
TNF- $\alpha$	628	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		14 JUN 2021