



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0294-2

**Lot Number:** A0080317

**Expiration Date:** 31 MAR 2027

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	138	pg/mL
IL-1 $\beta$	1990	pg/mL
IL-2	314	pg/mL
IL-4	775	pg/mL
IL-5	337	pg/mL
IL-6	51500	pg/mL
KC/GRO	803	pg/mL
IL-10	2720	pg/mL
IL-13	1600	pg/mL
TNF- $\alpha$	1080	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		22 APR 2022