

**CERTIFICATE OF ANALYSIS**


**Product Description:** Proinflammatory Panel 1 (rat) Calibrator Blend  
**Catalog Number:** C0294-2  
**Lot Number:** A0080410  
**Expiration Date:** 30 JUN 2030  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	438	pg/mL
IL-1 $\beta$	1530	pg/mL
IL-2	761	pg/mL
IL-4	535	pg/mL
IL-5	453	pg/mL
IL-6	94500	pg/mL
KC/GRO	997	pg/mL
IL-10	7150	pg/mL
IL-13	1460	pg/mL
TNF- $\alpha$	1210	pg/mL

**Reconstitution Volume:** 1000  $\mu$ L

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. 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 AUG 2025