

**CERTIFICATE OF ANALYSIS**


**Product Description:** Proinflammatory Panel 1 (rat) Calibrator Blend  
**Catalog Number:** C0294-2  
**Lot Number:** A0080380  
**Expiration Date:** 31 MAR 2029  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	456	pg/mL
IL-1 $\beta$	2180	pg/mL
IL-2	758	pg/mL
IL-4	544	pg/mL
IL-5	478	pg/mL
IL-6	84500	pg/mL
KC/GRO	1110	pg/mL
IL-10	6610	pg/mL
IL-13	1540	pg/mL
TNF- $\alpha$	1260	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		08 MAY 2024