



### CERTIFICATE OF ANALYSIS

**Product Description:** Proinflammatory Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0049-2  
**Lot Number:** A0080385  
**Expiration Date:** 31 DEC 2028  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	1380	pg/mL
IL-1 $\beta$	615	pg/mL
IL-2	1380	pg/mL
IL-4	181	pg/mL
IL-6	823	pg/mL
IL-8	583	pg/mL
IL-10	376	pg/mL
IL-12p70	457	pg/mL
IL-13	503	pg/mL
TNF- $\alpha$	354	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		28 FEB 2024