



### CERTIFICATE OF ANALYSIS

**Product Description:** Proinflammatory Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0049-2  
**Lot Number:** A0080368  
**Expiration Date:** 30 JUN 2028  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	1450	pg/mL
IL-1 $\beta$	642	pg/mL
IL-2	1530	pg/mL
IL-4	238	pg/mL
IL-6	728	pg/mL
IL-8	622	pg/mL
IL-10	376	pg/mL
IL-12p70	500	pg/mL
IL-13	560	pg/mL
TNF- $\alpha$	393	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		25 JUL 2023