

**CERTIFICATE OF ANALYSIS**


**Product Description:** Proinflammatory Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0049-2  
**Lot Number:** A0080442  
**Expiration Date:** 31 AUG 2030  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	1600	pg/mL
IL-1 $\beta$	562	pg/mL
IL-2	1510	pg/mL
IL-4	245	pg/mL
IL-6	730	pg/mL
IL-8	557	pg/mL
IL-10	348	pg/mL
IL-12p70	457	pg/mL
IL-13	517	pg/mL
TNF- $\alpha$	346	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		02 OCT 2025