

**CERTIFICATE OF ANALYSIS**


**Product Description:** Proinflammatory Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0049-2  
**Lot Number:** A0080428  
**Expiration Date:** 30 APR 2030  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	1490	pg/mL
IL-1 $\beta$	605	pg/mL
IL-2	1600	pg/mL
IL-4	258	pg/mL
IL-6	809	pg/mL
IL-8	571	pg/mL
IL-10	404	pg/mL
IL-12p70	507	pg/mL
IL-13	583	pg/mL
TNF- $\alpha$	356	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		29 MAY 2025