

**CERTIFICATE OF ANALYSIS**


**Product Description:** Proinflammatory Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0049-2  
**Lot Number:** A0080433  
**Expiration Date:** 30 APR 2030  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	1190	pg/mL
IL-1 $\beta$	479	pg/mL
IL-2	1580	pg/mL
IL-4	277	pg/mL
IL-6	764	pg/mL
IL-8	548	pg/mL
IL-10	360	pg/mL
IL-12p70	509	pg/mL
IL-13	592	pg/mL
TNF- $\alpha$	361	pg/mL

**Reconstitution Volume:** 1000  $\mu$ L

This product was manufactured and tested according to MSD documents. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Michelle Likiam	Quality		19 JUN 2025