

**CERTIFICATE OF ANALYSIS**

**Product Description:** Proinflammatory Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0049-2  
**Lot Number:** A0080378  
**Expiration Date:** 31 AUG 2028  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	1370	pg/mL
IL-1 $\beta$	627	pg/mL
IL-2	1510	pg/mL
IL-4	230	pg/mL
IL-6	712	pg/mL
IL-8	588	pg/mL
IL-10	377	pg/mL
IL-12p70	443	pg/mL
IL-13	545	pg/mL
TNF- $\alpha$	376	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		03 OCT 2023