



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080407
Expiration Date: 31 AUG 2029
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1350	pg/mL
IL-1 β	546	pg/mL
IL-2	1250	pg/mL
IL-4	195	pg/mL
IL-6	809	pg/mL
IL-8	591	pg/mL
IL-10	375	pg/mL
IL-12p70	481	pg/mL
IL-13	527	pg/mL
TNF- α	360	pg/mL

Reconstitution Volume: 1000 μ 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
Review/Approval	Karen Hamilla	Quality		10 OCT 2024