



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080363
Expiration Date: 31 MAR 2028
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1360	pg/mL
IL-1 β	606	pg/mL
IL-2	1500	pg/mL
IL-4	242	pg/mL
IL-6	721	pg/mL
IL-8	679	pg/mL
IL-10	374	pg/mL
IL-12p70	472	pg/mL
IL-13	577	pg/mL
TNF- α	368	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		25 APR 2023