



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080365
Expiration Date: 30 APR 2028
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1520	pg/mL
IL-1 β	650	pg/mL
IL-2	1580	pg/mL
IL-4	269	pg/mL
IL-6	725	pg/mL
IL-8	593	pg/mL
IL-10	393	pg/mL
IL-12p70	500	pg/mL
IL-13	637	pg/mL
TNF- α	375	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		02 JUN 2023