



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080382
Expiration Date: 30 NOV 2028
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1340	pg/mL
IL-1 β	625	pg/mL
IL-2	1440	pg/mL
IL-4	218	pg/mL
IL-6	822	pg/mL
IL-8	586	pg/mL
IL-10	365	pg/mL
IL-12p70	478	pg/mL
IL-13	551	pg/mL
TNF- α	369	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		20 DEC 2023