



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080397
Expiration Date: 31 MAY 2029
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1400	pg/mL
IL-1 β	630	pg/mL
IL-2	1470	pg/mL
IL-4	220	pg/mL
IL-6	816	pg/mL
IL-8	575	pg/mL
IL-10	382	pg/mL
IL-12p70	497	pg/mL
IL-13	514	pg/mL
TNF- α	362	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		14 JUN 2024