



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080395
Expiration Date: 31 MAR 2029
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1330	pg/mL
IL-1 β	632	pg/mL
IL-2	1450	pg/mL
IL-4	215	pg/mL
IL-6	820	pg/mL
IL-8	595	pg/mL
IL-10	378	pg/mL
IL-12p70	476	pg/mL
IL-13	531	pg/mL
TNF- α	363	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		13 MAY 2024