



CERTIFICATE OF ANALYSIS

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend
Catalog Number: C0049-2
Lot Number: A0080400
Expiration Date: 30 JUN 2029
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	1060	pg/mL
IL-1 β	624	pg/mL
IL-2	1430	pg/mL
IL-4	183	pg/mL
IL-6	784	pg/mL
IL-8	575	pg/mL
IL-10	378	pg/mL
IL-12p70	442	pg/mL
IL-13	514	pg/mL
TNF- α	344	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		12 AUG 2024